

EARLY ACCESS PROTOCOL

Monitoring HiBiT-KRAS(WT) Levels in HiBiT-KRAS(WT) KI HEK293 Cells (CPM)

Instructions for Use of
CS3023186R

These materials provide early access to innovative solutions before a product is available in our catalog. They are sold "as is" with no warranty either stated or implied. Changes may be incorporated without notice. These materials are provided with a data report; however, they do not have other supporting documentation, such as a Certificate of Analysis.

Early Access materials are intended For Research Use Only.

Monitoring HiBiT-KRAS(WT) Levels in HiBiT-KRAS(WT) KI HEK293 Cells (CPM)

All technical literature is available at: www.promega.com/protocols/
Visit the website to verify that you are using the most current version of this Early Access Protocol.
Email Promega Technical Services if you have questions on use of this system: techserv@promega.com

1. Description	2
2. Product Components and Storage Conditions	2
3. Before You Begin	2
3.A. Media for Cell Propagation and Freezing	2
3.B. Cell Propagation	3
4. Nano-Glo® HiBiT Lytic Detection Assay Protocol	4
4.A. Preparing Cells for Nano-Glo® HiBiT Lytic Assay	4
4.B. Nano-Glo® HiBiT Lytic Detection Assay	4
4.C. CellTiter-Glo® 2.0 Cell Viability Assay	4
4.D. Calculating Signal-to-Background Ratios	5



1. Description

This early access material is designed for monitoring HiBiT-KRAS(WT) levels in HEK293 cells that have been genetically modified via CRISPR-Cas9 to fuse HiBiT to the N terminus of KRAS(WT). The resultant clone cells have been verified to be heterozygous for the HiBiT insertion.

2. Product Components and Storage Conditions

PRODUCT	SIZE	PART NUMBER
HiBiT-KRAS(WT) KI HEK293 Cells (CPM)	1 each	CS3023186R

Includes:

- 2 vials HiBiT-KRAS(WT) KI HEK293 Cells (CPM) (CS3023185)

Storage Conditions: Store the HiBiT-KRAS(WT) KI HEK293 Cells (CPM)^(a-e) at -140°C in the liquid nitrogen vapor phase, upon receipt.

3. Before You Begin

Materials to Be Supplied By the User

- white, 96-well plate (e.g., Costar Cat.# 3917)
- miscellaneous tissue culture reagents and equipment
- DPBS (1X) (e.g., GIBCO® Cat.# 14190-144)
- 0.25% Trypsin-EDTA (GIBCO® Cat.# 25200-056)
- CellTiter-Glo® 2.0 Cell Viability Assay (Cat.# G9241)
- Nano-Glo® HiBiT Lytic Detection System (Cat.# N3030, N3040, N3050), which includes: Nano-Glo® HiBiT Lytic Substrate; Nano-Glo® HiBiT Lytic Buffer; and LgBiT Protein
- DMEM (e.g., Thermo Fisher Scientific Cat.# 11995065)
- FBS (e.g., VWR Cat.# 89510-194)
- DMSO (e.g., Sigma Cat.# D2650)

3.A. Media for Cell Propagation and Freezing

cell propagation and assay medium

- 90% DMEM
- 10% FBS

freezing medium

- 85% DMEM
- 10% FBS
- 5% DMSO

3.B. Cell Propagation

Thawing Cells for Culture

1. Remove vial from long-term storage (-140°C or below) and incubate in 37°C water bath for 2–3 minutes until thawed.
2. Transfer cells (1ml) into a 15ml centrifuge tube containing 9ml of propagation medium prewarmed to 37°C .
3. Centrifuge for 5 minutes at $200 \times g$.
4. Remove medium and suspend pellet in 10ml of prewarmed propagation medium. Place cell suspension in a T75 tissue culture flask.
5. Culture in a humidified 37°C , 5% CO_2 incubator.
6. Maintain cells in propagation medium and passage when approximately 80% confluent. A subcultivation ratio of 1:3–1:10 is recommended.

Freezing/Banking Cells

7. Remove propagation medium and wash cells once with DPBS.
8. Remove DPBS and add an appropriate volume of 0.25% Trypsin-EDTA, such that the entire surface area of the tissue culture dish is covered.
9. Incubate at 37°C for 5 minutes or until cells can be easily removed from the surface of the tissue culture plate.
10. Add an equal volume of propagation medium to the tissue culture plate and transfer the cell suspension into a conical tube for centrifugation.
11. Centrifuge cells 5 minutes at $200 \times g$ and carefully remove supernatant.
12. Gently suspend cell pellet in cold freezing medium at 2×10^6 – 5×10^6 cells/ml and aliquot into appropriately-sized cryovials.
13. Freeze in a controlled-rate freezer.
14. Store long term in liquid nitrogen vapor phase.

4. Nano-Glo® HiBiT Lytic Detection Assay Protocol

See *Nano-Glo® HiBiT Lytic Detection System Technical Manual #TM516* and *CellTiter-Glo® 2.0 Assay Technical Manual #TM403* for detailed instructions.

4.A. Preparing Cells for Nano-Glo® HiBiT Lytic Assay

1. Adjust the cell density to 2×10^5 cells/ml in assay medium. The cell number can be adjusted to increase luminescent signal.
2. Dispense 100µl per well of cell suspension (2×10^4 cells) into a white 96-well plate.
3. Plate replicates for both the lytic assay and the cell viability assay at the same time. Do not place the lytic and viability samples in adjacent wells, as the luminescence from the CellTiter-Glo® 2.0 Reagent may be bright enough to affect readings in neighboring wells. If placing lytic and cell viability samples in adjacent wells is unavoidable, perform the Nano-Glo® HiBiT Assay first, read the luminescence, then perform the CellTiter-Glo® Assay.
4. Cells can be assayed the day of plating or later, depending on the type of plate and cell density.

4.B. Nano-Glo® HiBiT Lytic Detection Assay

Immediately prior to Nano-Glo® HiBiT Lytic measurements, prepare 2X Nano-Glo® HiBiT Lytic Detection Reagent by diluting LgBiT Protein 1:100 and Nano-Glo® HiBiT Lytic Substrate 1:50 into an appropriate volume of room temperature Nano-Glo® HiBiT Lytic Buffer. Prepare enough 2X detection reagent for the number of wells to be assayed, including extra volume to account for pipetting error (for example, number of wells + 10%).

1. Add 100µl per well of 2X Nano-Glo® HiBiT Lytic Detection Reagent for a 96-well plate. Mix plate on an orbital shaker for 10–20 minutes at 500rpm.
2. Measure luminescence on a GloMax® Discover System (Cat.# GM3000) or other plate-reading luminometer.

4.C. CellTiter-Glo® 2.0 Cell Viability Assay

1. Equilibrate CellTiter-Glo® 2.0 Reagent to room temperature prior to starting the assay.
2. During the Nano-Glo® HiBiT Lytic Assay, plate replicate wells to examine cell viability with a final volume of 100µl per well.
3. Add 100µl per well of CellTiter-Glo® 2.0 Reagent for a 96-well plate. Mix plate on an orbital shaker for 10–20 seconds at 350rpm.
4. Measure luminescence on a GloMax® Discover System or other luminometer.

4.D. Calculating Signal-to-Background Ratios

$$\text{Signal-to-background ratio} = \frac{S/\text{CTG, S}}{P/\text{CTG, P}}$$

where:

S = sample reporter signal

P = parental/control cells reporter signal

CTG, S = sample CellTiter-Glo® 2.0 signal

CTG, P = parental/control cells CellTiter-Glo® 2.0 signal

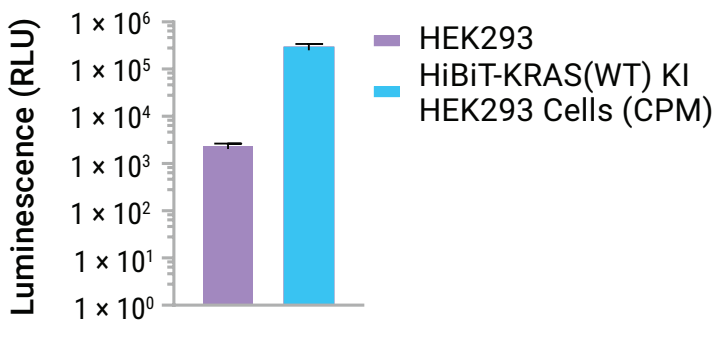


Figure 1. Measurement of endogenously-tagged HiBiT-KRAS(WT) by the Nano-Glo® HiBiT Lytic Detection Assay.

HEK293 parental cells and HiBiT-KRAS(WT) KI HEK293 Cells (CPM) were plated into a white 96-well assay plate at 20,000 cells/well in triplicate. HiBiT-KRAS(WT) levels were measured by adding an equal volume of 2X Nano-Glo® Lytic Assay Detection Reagent. Luminescence was measured using a GloMax® Discover System. Relative light units (RLU) are shown on the y axis. Normalizing to CellTiter-Glo®, the signal-to-background ratio is approximately 156.6 after thawing (data not shown). Data were graphed and analyzed using GraphPad Prism® software.



^(a)Limited Use Label License

BY USE OF THIS MATERIAL, RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE LABEL LICENSE.

Researchers may use this material for research use only. No transfer or commercial use of this material, progeny of this material, or derivatives of this material is allowed. Commercial use means any and all uses of this material or derivatives by a party in exchange for consideration, including, but not limited to (1) use in further product manufacture; (2) use in provision of services, information or data; and (3) resale of the material or its derivatives, whether or not such material or derivatives are resold for use in research.

Researchers may modify the material, except researchers shall have no right to modify or otherwise create variations of the nucleotide sequences of the subunits of NanoBiT[®] Technology, except that researchers may create fused gene sequences at the termini of the subunits of NanoBiT[®] Technology. No other use of this material or its derivatives is authorized without the prior express written consent of Promega.

In addition, researchers must either:

(4a) use Nano-Glo[®]-branded luminescent assay reagents (LARs) for all determinations of luminescence activity of this material and its derivatives, or

(4b) contact Promega to obtain a license for use of the material and its derivatives with LARs not manufactured by Promega.

For uses of Nano-Glo[®]-branded LARs intended for energy transfer (such as bioluminescence resonance energy transfer) to acceptors other than a genetically encoded autofluorescent protein, researchers must either:

(5a) use NanoBRET[™]-branded energy acceptors (e.g., BRET-optimized HaloTag[®] ligands) for all determinations of energy transfer activity by this material and its derivatives, or

(5b) contact Promega to obtain a license for use of the material and its derivatives for energy transfer assays to energy acceptors not manufactured by Promega.

With respect to any uses outside this label license, including any diagnostic, therapeutic, prophylactic or commercial uses, please contact Promega for supply and licensing information. PROMEGA MAKES NO REPRESENTATIONS OR WARRANTIES OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING FOR MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, WITH REGARD TO THE MATERIAL. The terms of this label license shall be governed under the laws of the State of Wisconsin, USA.

^(b)Broad Limited License

By purchase of this material, Promega Corporation ("Promega") conveys to purchaser (the "Limited Licensee") the non-transferable right to use the material purchased from Promega and derivatives made by Limited Licensee, solely for research conducted by Limited Licensee in accordance with the following

requirements: (i) Limited Licensee shall not sell or otherwise transfer material or derivatives to any other person or entity, or use material or derivatives to perform services for the benefit of any other person or entity, except to transfer the material or derivatives to a Limited Licensee affiliate provided such affiliate agrees to be bound by the terms of this Limited License (ii) Limited Licensee shall use material and derivatives only for its internal research within the Customer Field, which may include internal research within the Customer Field in connection with product research, and not for any Commercial Purposes, (iii) Limited Licensee shall use material and derivatives in compliance with all applicable laws and regulations, including without limitation applicable human health and animal welfare laws and regulations, (iv) Institutions shall provide no warranties of any kind to Limited Licensee (statutory or implied) concerning the Intellectual Property or material or derivatives, including without limitation, as to product quality, condition, description, merchantability, fitness for a particular purpose, noninfringement of intellectual property rights or the absence of latent or other defects, and all such warranties are hereby expressly disclaimed, (v) Institutions shall expressly disclaim any warranty regarding results obtained through the use of material or derivatives, including without limitation any claim of inaccurate, invalid or incomplete results, (vi) Institutions and their directors, trustees, officers, employees, agents, faculty, affiliated investigators, and students, shall have no liability to Limited Licensee, including, without limitation, for any loss of use or profits, business interruption or any consequential, incidental, special or other indirect damages of any kind, regardless of how caused and regardless of whether an action in contract, tort, strict product liability or otherwise, (vii) Limited Licensee shall indemnify, defend and hold harmless the Indemnitees and HHMI Indemnitees against any liability, damage, loss, or expense (including without limitation reasonable attorneys' fees and expenses) incurred by or imposed upon any of the Indemnitees or HHMI Indemnitees, as applicable, in connection with any claims, suits, investigations, actions, demands or judgments arising out of or related to the exercise of any rights granted to Limited Licensee under this limited license or any breach of limited license by such Limited Licensee, provided that, to the extent the foregoing is not permitted by law, Limited Licensee agrees, to the extent permitted by law, that it, and not the Indemnitees or HHMI Indemnitees, as applicable, shall be responsible for any liability, damage, loss or expense arising out of or related to the exercise of any rights granted to Limited Licensee under this limited license or any breach of this limited license by Limited Licensee, and (viii) material and derivatives and their use may be the subject of one or more issued patents or pending patent applications owned by one or more Institutions and the purchase of material does not convey a license under any claims in the foregoing patents or patent applications directed to the material or derivatives or use, production or commercialization thereof, except as expressly set forth in this limited license. Nothing in this limited license shall be construed to confer any rights upon Limited Licensee by implication, estoppel, or otherwise as to any technology or patent rights of Broad or any other entity other than the Intellectual Property. In addition, nothing in this limited license shall be construed to confer upon Limited Licensee or any third party any rights under or to the Intellectual Property outside of the Customer Field.

Commercial Purposes means (a) the practice, performance or provision of any methods, process or service, or (b) the manufacture, production, sale, use, distribution, disposition or importing of any product, in each case (a) or (b) for consideration of any kind, for the purpose of sale or commercial exploitation, or on any other commercial basis.

Customer Field means use as a research tool for research purposes and use for quality control/quality assurance testing and product lot release assays; provided, however, that notwithstanding the foregoing, the Customer Field shall expressly exclude:

- 1) Any human or clinical use, including, without limitation, any administration into humans or any diagnostic or prognostic use;
- 2) Any human germline modification, including modifying the DNA of human embryos or human reproductive cells;
- 3) Any in vivo veterinary or livestock use;
- 4) The development, manufacture (including any bioproduction), distribution, importation, exportation, transportation, sale, offer for sale, marketing, promotion or other exploitation or use of the Intellectual Property or the material or derivatives as a therapeutic or diagnostic for humans or animals;
- 5) Products that provide nutritional benefits and are regulated by a regulatory authority as a drug or biologic pursuant to Section 505 of the Federal Food, Drug, and Cosmetic Act of 1938, as amended, Section 351 of the Public Health Service Act of 1944, as amended, or any successor laws, or equivalent laws or regulations in jurisdiction outside the United States;
- 6) Any agricultural use, including but not limited to, the use or application in the cultivation, growth, manufacture, exportation, or production of any tobacco product; and
- 7) Any use or application relating to gene drive.

HHMI Indemnities means Howard Hughes Medical Institute and its trustees, officers, employees, and agents.

Indemnitees means each Institution, their affiliates, and their current and former trustees, directors, officers, faculty, affiliated investigators, students, employees, medical and professional staff and agents and their respective successors, heirs and assigns.

Institutions means each of The Broad Institute, Inc., Massachusetts Institute of Technology, Harvard College, University of Iowa Research Foundation, University of Tokyo, The Rockefeller University, New York University, New York Genome Center, and Whitehead Institute of Biomedical Research collectively.

Intellectual Property means (a) any patents and patent applications listed on Appendix A; (b) any substitutions, divisions, continuations, reissues, renewals, re-examinations or extensions of any patents or patent applications of (a), any continuations-in-part of any patents or patent applications of (a) to the extent the claims are directed to subject matter specifically described in a patent or patent application set forth on Appendix A, and any application claiming priority to a provisional application set forth on Appendix A to the extent the claims are directed to subject matter specifically described in a provisional patent application set forth on Appendix A; (c) any letters patents and/or the equivalent issuing therefrom in any jurisdiction based on, and only to the extent of, any of the foregoing; and (d) any foreign or international equivalents of any of the foregoing, including without limitation any applications filed under the Patent Cooperation Treaty.



©Sigma-Aldrich CRISPR Use License Agreement

This Product and its use are the subject of one or more of the following patents and applications controlled by Sigma-Aldrich Co., LLC (SIGMA): Patent applications and issued patents that entered their respective National Stage from PCT International Pub. No. WO 2014/089290, including, but not limited to, the following, and substitutions, divisions, continuations, continuations-in-part, reissues, reexaminations, and extensions thereof: Australia Patent Nos. 2013355214; 2017204031; and 2018229489; Canada Patent No. 2,891,347; China Patent No. CN105142669; European Patent Nos. EP 3 138 910 B1, 3 138 911 B1, and EP 3 138 910 B1; Israel Patent No. IL238856; Singapore Patent No. 11201503824S; South Korea Patent No. 10-1844123; and U.S. Patent Application Serial Nos. 15/188,911; 15/188,924; 15/188,927; 15/188,931; and 15/456,204 (the "Patent Rights").

BEFORE OPENING OR USING THIS PRODUCT, PLEASE READ THE TERMS AND CONDITIONS SET FORTH IN THIS LICENSE AGREEMENT. YOUR USE OF THIS PRODUCT SHALL CONSTITUTE ACKNOWLEDGMENT AND ACCEPTANCE OF THESE TERMS AND CONDITIONS. If you do not agree to use this Product pursuant to the terms and conditions set out in this License Agreement, please contact Promega Corporation within ten (10) days of receipt to return the unused and unopened Product for a full refund; provided, however, that custom-made Products may not be returned for a refund.

The purchase of this Product conveys to you (the "Buyer") the NON-TRANSFERABLE right to use the Product for Licensed Research Use (see definition below) subject to the conditions set out in this License Agreement. If you wish to use this Product for any purpose other than Licensed Research Use, you must first obtain an appropriate license (see information set out below).

This Product may not be used for any purpose other than Licensed Research Use. Your right to use this Product for Licensed Research Use is subject to the following conditions and restrictions:

1. "Licensed Research Use" means any use for research purposes, with the following exceptions: (i) Buyer may not sell or otherwise transfer the Product (including without limitation any material that contains the Product in whole or part) or any Related Material to any other third party; (ii) Buyer may use the Product, components of the Product, and any Related Material, only for your internal research within the Field, and not for any Commercial Purposes (for any use of Product, components of a Product and any Related Material by any entity, including but not limited to contract service providers ("CRO") offering services for consideration, please contact Promega for additional information) and this License Agreement does not grant or in any way convey any rights to Buyer to further edit Products, components of Products or Related Materials under any claims in the foregoing patents and patent applications; (iii) Buyer shall use the Product and any Related Material in compliance with all applicable laws and regulations, including without limitation applicable human health and animal welfare laws and regulations; (iv) SIGMA shall provide no warranties of any kind to the Buyer (statutory or implied) concerning the Patent Rights, the Product or any Related Materials, including without limitation, as to product quality, condition, description, merchantability, fitness for a particular purpose, noninfringement of intellectual property rights or the absence of latent or other defects, and all such warranties are hereby expressly disclaimed; (v) SIGMA shall expressly disclaim any warranty regarding results obtained through the use of the Product or Related Materials, including without limitation any claim of inaccurate, invalid or incomplete results; (vi) SIGMA, and their directors, officers, employees and agents, shall have no liability to the Buyer, including, without limitation, for any loss of use or profits, business interruption or any consequential, incidental, special or other indirect damages of any kind, regardless of how caused and regardless of whether an action in contract, tort, strict product liability or otherwise; (vii) the Buyer shall indemnify, defend, and hold harmless SIGMA and their current and former directors, officers, employees and agents, and their respective successors, heirs and assigns (the "Indemnitees") against any liability, damage, loss, or expense (including without limitation reasonable attorneys' fees and expenses) incurred by or imposed upon any of the Indemnitees in connection with any claims, suits, investigations, actions, demands or judgments arising out of or related to the exercise of any rights granted to the Buyer hereunder or any breach of this License Agreement by such Buyer; provided that, to the extent the foregoing is not permitted by law, the Buyer agrees, to the extent permitted by law, that it, and not the Indemnitees shall be responsible for any liability, damage, loss or expense arising out of or related to the exercise of any rights granted to the Buyer under this License Agreement or any breach of this License Agreement by Buyer, and (viii) the Product and its creation and/or use may be the subject of one or more issued patents and/or pending patent applications owned by SIGMA and the purchase of the Product and its use as defined in this License Agreement does not convey to the Buyer any license under any claims in the foregoing patents or patent applications except for the limited rights as described in this License Agreement directed toward use of the Licensed Product or commercialization thereof.

2. For purposes of Section 1 above, the following definitions shall apply:

"Commercial Purposes" means (a) the practice, performance or provision of any method, process or service or (b) the manufacture, sale, distribution, disposition or importing of any product, in each case (a) or (b) for consideration, e.g., a fee, or on any other commercial basis.

"Field" means use as a research tool for research purposes which includes, but is not limited to drug discovery and development, quality control/quality assurance testing and product lot release assays; provided, however, that notwithstanding the foregoing, the Field shall expressly exclude (a) any in vivo and ex vivo human or clinical use, including, without limitation, any administration into humans or any diagnostic or prognostic use, (b) the creation of transgenic rodent models and/or derivatives thereof (such derivatives including, but not limited to, rodents' cells and rodents' organs), (c) any in vivo veterinary or livestock use, or non-research agricultural use, or (d) the manufacture, distribution, importation, exportation, transportation, sale, offer for sale, marketing, promotion or other exploitation or use of, or as, a therapeutic or diagnostic for humans or animals.

"Related Materials" means any progeny, modification or derivative of a Product.

3. Your right to use the Product will terminate immediately if you fail to comply with these terms and conditions.

For information on purchasing a license a) to this Product for purposes other than Licensed Research Use, or b) to the foregoing patents and patent applications generally, contact your local SIGMA sales representative, who will refer you to the proper licensing representative, or in the USA call 800-325-3010.

^(d)Caribou Biosciences Limited Use Label License

BEFORE OPENING OR USING THIS PRODUCT, PLEASE READ THE TERMS AND CONDITIONS SET FORTH IN THIS LIMITED USE LABEL LICENSE. YOUR USE OF THIS PRODUCT SHALL CONSTITUTE ACKNOWLEDGMENT AND ACCEPTANCE OF THESE TERMS AND CONDITIONS. If you do not agree to use this Product pursuant to the terms and conditions set out in this Limited Use Label License, please contact Promega Corporation within ten (10) calendar days of receipt to return the unused and unopened Product for a full refund; provided, however, that custom-made Products may not be returned for a refund.

Promega is selling this Product under a research use only license from Caribou Biosciences, Inc. under certain patents and patent applications owned or controlled by Caribou. For information on obtaining rights beyond those set forth in this Limited Use Label License please contact licensing@promega.com or Promega Corporation, 2800 Woods Hollow Road, Madison, WI 53711 USA, Attn: Licensing.

The purchase of this Product conveys to you, the Buyer, the NON-TRANSFERABLE right to use this Product solely for Research Use, subject to the conditions set out in this Limited Use Label License. If you wish to use this Product for any purpose other than Research Use, you must first obtain an appropriate license. This Product may not be used for any purpose other than Research Use. Your right to use this Product for Research Use is subject to the following conditions and restrictions:

1. Research Use means any use for internal research purposes, (e.g., for screening, characterization, quality control/quality assurance testing, product lot release assays, and/or toxicology purposes) for the sole benefit of you, the Buyer.

2. Research Use specifically excludes the following:

(a) Sale or otherwise transfer of this Product, including without limitation any material that contains this Product in whole or part or any progeny, modification or derivative of this Product (collectively, "Related Materials") to any other third party; and

(b) Commercial purposes and prohibited activities including use of this Product and Related Materials in (i) any sale, marketing, distribution, disposition, importation, or exportation, whether for a fee or any other commercial basis; (ii) any commercial activity including, but not limited to, any use in manufacturing or bioproduction; (iii) preclinical or clinical testing or other activity directed toward the submission of data to the U.S. Food and Drug Administration, or any other regulatory agency in any country or jurisdiction where the active agent in such studies comprises this Product; (iv) use in a product for, or for any use in, human or animal therapeutics, diagnostics, or prophylactics; (v) activity in an agricultural field trial or any activity directed toward the submission of data to the U.S. Department of Agriculture or any other agriculture regulatory agency; (vi) modification of the human germline, including editing human embryos or reproductive cells; (vii) stimulation of biased inheritance of a particular gene or trait or set of genes or traits; and/or (viii) use of self-administration.

3. You, the Buyer, shall use this Product and Related Materials in compliance with all applicable laws and regulations.

4. No warranties of any kind are provided to the you, the Buyer, (statutory or implied) concerning this Product or Related Materials, including without limitation, as to product quality, condition, description, merchantability, fitness for a particular purpose, noninfringement of intellectual property rights, or the absence of latent or other defects, and all such warranties are hereby expressly disclaimed.

5. Your right to use this Product and Related Materials will terminate immediately if you fail to comply with these terms and conditions.

^(e) U.S. Pat. Nos. 9,797,890 and 10,648,971 and other patents and patents pending.

© 2024 Promega Corporation. All Rights Reserved.

CellTiter-Glo, GloMax and Nano-Glo are registered trademarks of Promega Corporation.

GIBCO is a registered trademark of Life Technologies Corp. GraphPad Prism is a registered trademark of GraphPad Software, Inc.

Materials may be covered by pending or issued patents or may have certain limitations. Please visit our website for more information.

Pricing and availability are subject to change without prior notice.