



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet
GITR:GITRL TR-FRET Assay Kit
Catalog # 79054

DESCRIPTION:

The GITR:GITRL TR-FRET Assay is designed to measure the inhibition of GITR binding to GITRL in a homogeneous 384 reaction format. This FRET-based assay requires no time-consuming washing steps, making it especially suitable for high throughput screening applications. The assay procedure is straightforward and simple; a sample containing biotinylated GITR, GITRL, anti-His Tb donor, dye-labeled acceptor, and an inhibitor is incubated for two hours. Then, the fluorescence intensity is measured using a fluorescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
71256	GITR, biotinylated	15 µg	-80°C	(Avoid freeze/thaw cycles!)
71190	GITRL-His	2 µg	-80°C	
	Anti-His Tb Donor	2 x 10 µl	-20°C	
	Dye-labeled Acceptor	2 x 10 µl	-20°C	
	3x GITR TR FRET Assay Buffer	4 ml	-20°C	
	White, non-binding, low volume, 384-well microtiter plate	1	Room temp.	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescence microplate reader capable of measuring Time Resolved Fluorescence Resonance Energy Transfer (TR-FRET)
Adjustable micropipettor and sterile tips

APPLICATIONS: Great for screening small molecular inhibitors for drug discovery and HTS applications.

STABILITY: At least 6 months from date of receipt when stored as directed.

REFERENCES:

1. Lechner, M.G., *et al. Immunotherapy*. 2011 Nov; **3(11)**: 1317–1340.
2. Ko, K., *et al., J. Exp. Med.* 2005; **202(7)**: 885-891.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

All samples and controls should be tested in at least duplicate.

Protocol for GITR assay

- 1) Dilute one part **3x GITR TR-FRET Assay Buffer** with 2 parts distilled water (3-fold dilution) to make **1x GITR Assay Buffer**. Make only a sufficient quantity needed for the assay; store remaining stock solution in aliquots at -20°C.
- 2) Dilute **Anti-His Tb Donor** 100-fold in **1x GITR Assay Buffer**. Make only sufficient quantities needed for the assay; store remaining stock solution in aliquots at -20°C.
- 3) Dilute **Dye-labeled acceptor** 100-fold in **1x GITR Assay Buffer**. Make only sufficient quantities needed for the assay; store remaining stock solution in aliquots at -20°C.
- 4) Thaw **GITR, Biotinylated** on ice. Upon first thaw, briefly spin tube containing **GITR, Biotinylated** to recover the full contents of the tube. Aliquot into single-use aliquots. Store remaining undiluted **GITR** at -80°C immediately. *Note: **GITR, Biotinylated** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 5) Dilute **GITR, Biotinylated** in **1x GITR Assay Buffer** to 8 µg/ml. Make only sufficient quantities needed for the assay; store remaining stock solution in aliquots at -20°C.
- 6) Prepare the master mixture: N wells x (3 µl diluted **GITR, Biotinylated** + 5 µl diluted **Anti-His Tb Donor** + 5 µl diluted **Dye-labeled acceptor**). Add 13 µl to every well.
- 6) Add 2 µl of inhibitor solution to each well designated "Test Inhibitor." Add 2 µl of the same solution without inhibitor (inhibitor buffer) to the wells labeled "Negative Control" and "Positive Control."
- 7) Add 5 µl **1x GITR Assay Buffer** to wells designated for "Negative Control."

	Positive Control	Negative Control	Test Inhibitor
GITR, Biotinylated (diluted)	3 µl	3 µl	3 µl
Anti-His Tb Donor	5 µl	5 µl	5 µl
Dye-labeled Acceptor	5 µl	5 µl	5 µl
Test Inhibitor	-	-	2 µl
Inhibitor Buffer (no inhibitor)	2 µl	2 µl	-
1x GITR Assay Buffer	-	5 µl	-
GITRL-His (diluted)	5 µl	-	5 µl
Total	20 µl	20 µl	20 µl

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 8) Thaw **GITRL-His** protein on ice. Upon first thaw, briefly spin tube containing protein to recover the full contents of the tube. Aliquot **GITRL-His** into single-use aliquots. Store remaining undiluted **GITRL-His** in aliquots at -80°C immediately. *Note: **GITRL-His** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 9) Dilute **GITRL-His** in **1x GTR Assay Buffer** to $0.2\ \mu\text{g/ml}$. Initiate reaction by adding $5\ \mu\text{l}$ of diluted **GITRL-His** to wells designated for the "Positive Control" and "Test Inhibitor." Discard any remaining diluted **GITRL-His** protein after use.
- 10) Incubate at room temperature for 1.5 hours.
- 11) Read the fluorescent intensity in a microtiter-plate reader capable of TR-FRET.

Instrument Settings

Reading Mode	Time Resolved
Excitation Wavelength	$320\pm 10\ \text{nm}$
Emission Wavelength	$620\pm 10\ \text{nm}$
Lag Time	$60\ \mu\text{s}$
Integration Time	$500\ \mu\text{s}$
Excitation Wavelength	$320\pm 20\ \text{nm}$
Emission Wavelength	$665\pm 10\ \text{nm}$
Lag Time	$60\ \mu\text{s}$
Integration Time	$500\ \mu\text{s}$

CALCULATING RESULTS:

Two sequential measurements should be conducted. Tb-donor emission should be measured at $620\ \text{nm}$ followed by dye-acceptor emission at $665\ \text{nm}$. Data analysis is performed using the TR-FRET ratio ($665\ \text{nm}$ emission/ $620\ \text{nm}$ emission).

If desired, data can be normalized to percent inhibition. Typically for inhibitor screens, the FRET value from the positive control is set to zero percent inhibition and the FRET value from the negative control is set to one hundred percent inhibition.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com

EXAMPLE OF ASSAY RESULTS:

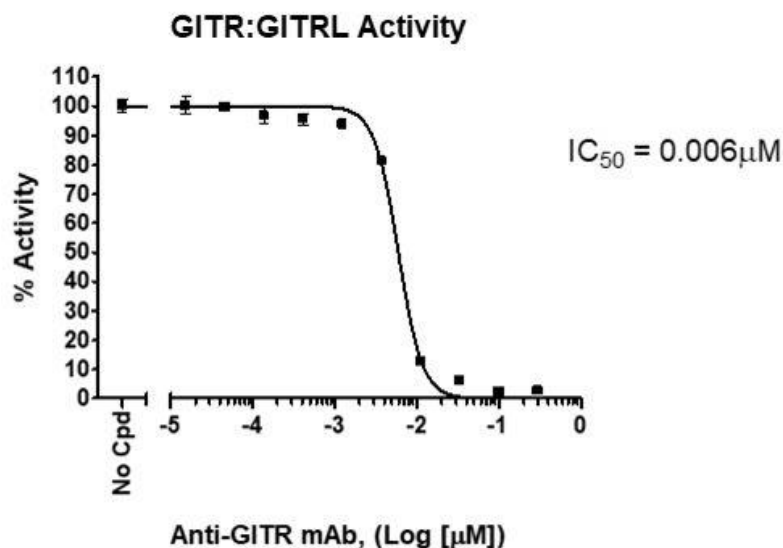


Figure Legend: Inhibition of GITR:GITRL interaction with anti-GITR antibody (BPS Bioscience, #79053). Data in the above graphs are expressed as FRET ratios. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com*

RELATED PRODUCTS:

<u>Product</u>	<u>Catalog #</u>	<u>Size</u>
GITR - HEK293 Recombinant Cell Line	79092	2 vials
GITR / NF-κB Luciferase Reporter (Luc) - Jurkat Cell Line	60546	2 vials
GITRL CHO-K1 Recombinant Cell Line	60547	2 vials
GITRL:GITR[Biotinylated] Inhibitor Screening Assay Kit	72061	96 rxns.
Anti-GITR Antibody	79053	100 µg
Anti-GITR Antibody, PE-labeled	71295-1	50 µg
Anti-GITR Antibody, PE-labeled	71295-2	100 µg
GITRL, His-tag (Human)	71190	100 µg
GITR (CD357), Fc Fusion, Biotin-labeled (Human)	71256	50 µg
GITR (CD357), Fc fusion (Human)	71172	100 µg

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com