



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

Data Sheet
RANK:RANKL TR-FRET Assay Kit
Catalog #79101-1
Size: 96 reactions

DESCRIPTION: The RANK:RANKL TR-FRET Assay is designed to measure the inhibition of RANK binding to RANKL. 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 RANK, RANKL, anti-His Tb donor, dye-labeled acceptor, and an inhibitor is incubated for one hour. Then, the fluorescence intensity is measured using a fluorescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
70822	RANK, Fc fusion (IgG1), Biotin Labeled (Human)	5 µg	-80°C	Avoid multiple freeze/thaw cycles!
71051	RANKL, His-Tag (Human)	2 µg	-80°C	
30017	Anti-His Tb Donor	3 x 10 µl	-20°C	
	Dye-labeled Acceptor	3 x 10 µl	-20°C	
79311	3x Immuno Buffer 1	4 ml	-20°C	
	White, 96-well half-area microtiter plate	1 unit	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: 12 months from date of receipt, when stored as recommended.

REFERENCES:

Annis, A., *et al. J. Amer. Chem. Soc.* 2004, **126(4)**: 15495-15503
Yan, T., *et al. J. Cell. Biochem.* 2001, **83(2)**: 320-325

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

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 W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Protocol for RANK assay

- 1) Dilute one part **3x Immuno Buffer 1** with 2 parts distilled water (3-fold dilution) to make **1x Immuno Buffer 1**. 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 Immuno Buffer 1**. 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 Immuno Buffer 1**. Make only sufficient quantities needed for the assay; store remaining stock solution in aliquots at -20°C.
- 4) Thaw **RANK, Biotinylated** on ice. Upon first thaw, briefly spin tube containing **RANK, Biotinylated** to recover the full contents of the tube. Aliquot into single-use aliquots. Store remaining undiluted **RANK** at -80°C immediately. Note: **RANK, Biotinylated** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.
- 5) Dilute **RANK, Biotinylated** in **1x Immuno Buffer 1** to 1.65 ng/μl. 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 (15 μl diluted **RANK, Biotinylated** + 25 μl diluted **Anti-His Tb Donor** + 25 μl diluted **Dye-labeled acceptor**). Add 65 μl to every well.
- 7) Add 10 μl of inhibitor solution to each well designated "Test Inhibitor." Add 10 μl of 10% DMSO in water inhibitor (inhibitor buffer) to the wells labeled "Negative Control" and "Positive Control."
- 8) Add 25 μl **1x Immuno Buffer 1** to wells designated for "Negative Control."

	Positive Control	Blank	Test Inhibitor
RANK, Biotinylated (1.65 ng/μl)	15 μl	15 μl	15 μl
Anti-His Tb Donor	25 μl	25 μl	25 μl
Dye-labeled Acceptor	25 μl	25 μl	25 μl
Test Inhibitor	-	-	10 μl
10% DMSO in water (Inhibitor buffer)	10 μl	10 μl	-
1x Immuno Buffer 1	-	25 μl	-
RANKL-His (0.5 ng/μl)	25 μl	-	25 μl
Total	100 μl	100 μl	100 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 W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 9) Thaw **RANKL-His** protein on ice. Upon first thaw, briefly spin tube containing protein to recover the full contents of the tube. Aliquot **RANKL-His** into single-use aliquots. Store remaining undiluted **RANKL-His** in aliquots at -80°C immediately. Note: **RANKL-His** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.
- 10) Dilute **RANKL-His** in **1x Immuno Buffer 1** to 0.5 ng/ μl . Initiate reaction by adding 25 μl of diluted **RANKL-His** to wells designated for the "Positive Control" and "Test Inhibitor." Discard any remaining diluted **RANKL-His** protein after use.
- 11) Incubate at room temperature for 1 hour.
- 12) Read the fluorescent intensity in a microtiter-plate reader capable of TR-FRET.

Instrument Settings

Reading Mode	Time Resolved
Excitation Wavelength	320 \pm 10 nm
Emission Wavelength	620 \pm 10 nm
Lag Time	60 μs
Integration Time	500 μs
Excitation Wavelength	320 \pm 20 nm
Emission Wavelength	665 \pm 10 nm
Lag Time	60 μs
Integration Time	500 μs

CALCULATING RESULTS:

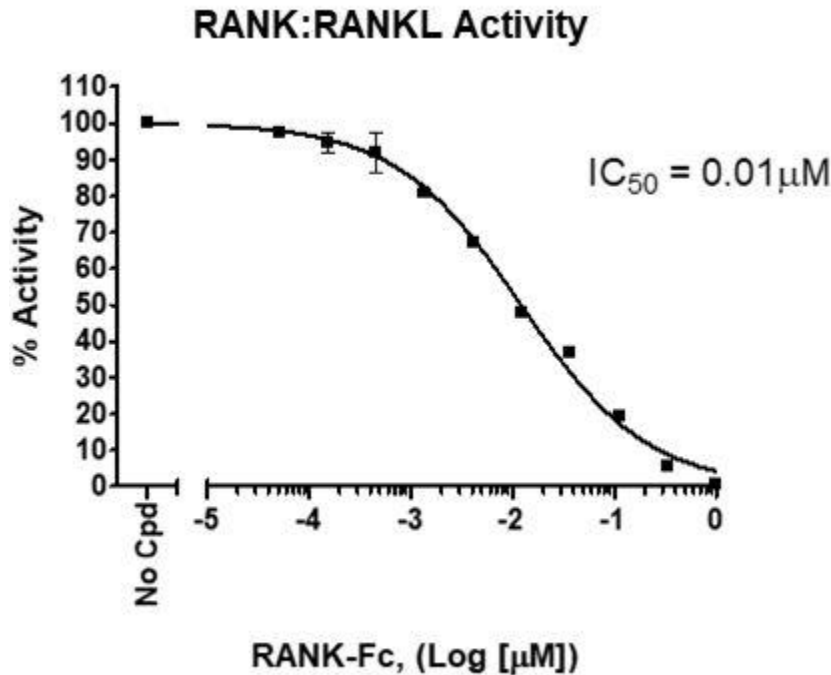
Two sequential measurements should be conducted. Tb-donor emission should be measured at 620 nm followed by dye-acceptor emission at 665 nm. Data analysis is performed using the TR-FRET ratio (665 nm emission/620 nm emission).

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:



Inhibition of RANK:RANKL interaction with unlabeled RANK (BPS Bioscience, #70823) and the RANK:RANKL TR-FRET Assay Kit (#79101-1). Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog#</u>	<u>Size</u>
RANKL, His-Tag (Human)	71051	100 µg
RANK, Fc fusion (IgG1), Avi-tag (Human)	70823	100 µg
RANK, Fc fusion (IgG1), Biotin Labeled (Human)	70822	50 µg
RANK:RANKL TR-FRET Assay	79101	384 reactions

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