



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: info@bpsbioscience.com

**Data Sheet**  
**L3MBTL1 TR-FRET Assay Kit**  
**Catalog # 55200**

**DESCRIPTION:** The L3MBTL1 TR-FRET Assay Kit is designed to measure the inhibition of L3MBTL1 binding to its substrate 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 terbium-labeled donor, dye-labeled acceptor, L3MBTL1, substrate, and an inhibitor is incubated for sixty minutes. Then, the fluorescence intensity is measured using a fluorescence reader.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
55000	GST-L3MBTL1(191-530)	10 µg	-80°C	<b>(Avoid freeze/ thaw cycles!)</b>
	Methylated MBT Ligand 1	50 µl	-80°C	
	Non-Methylated MBT Ligand 1	15 µl	-20°C	
	Tb-labeled donor	20 µl	-20°C	
	Dye-labeled acceptor	20 µl	-20°C	
	3x L3MBTL1 TR-FRET Assay Buffer	4 ml	-20°C	
Fisher 07-200-330	Nonbinding Corning, low volume, microtiter plate	1 plate	Room temp.	

**MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

Fluorescent 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.

**REFERENCE(S):** Min, J., *et al. Nat Struct Mol Biol* 2007; Dec;**14(12)**:1229-30.

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: info@bpsbioscience.com

## ASSAY PROTOCOL:

*All samples and controls should be tested in duplicate.*

### Protocol for L3MBTL1 assay

- 1) Dilute one part **3x L3MBTL1 TR-FRET Assay Buffer** with 2 parts distilled water (3-fold dilution) to make **1x L3MBTL1 Assay Buffer**. Make only a sufficient quantity needed for the assay; store remaining stock solution in aliquots at -20°C.
- 2) Dilute **Tb-labeled donor** and **Dye-labeled acceptor** 100-fold in **1x L3MBTL1 Assay Buffer**. Make only sufficient quantities needed for the assay; store remaining stock solution in aliquots at -20°C.
- 3) Add 5 µl of diluted **Tb-labeled donor**, and 5 µl of diluted **Dye-labeled acceptor** to each well designated "Test Inhibitor", "Substrate Control", and "Positive Control".
- 4) 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 "Substrate Control", and "Positive Control".

	<b>Substrate Control</b>	<b>Positive Control</b>	<b>Test Inhibitor</b>
Tb-labeled 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	–
Methylated MBT Ligand 1		5 µl	5 µl
Non-Methylated MBT Ligand 1	5 µl	–	–
L3MBTL1 3.5 ng/µl	3 µl	3 µl	3 µl
<b>Total</b>	<b>20 µl</b>	<b>20 µl</b>	<b>20 µl</b>

- 5) Thaw **Methylated MBT Ligand 1** and **Non-Methylated MBT Ligand 1** on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full contents of the tube. Aliquot each ligand into single-use aliquots. Store remaining undiluted ligand at -80°C immediately. *Note: each ligand is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots.*
- 6) Individually dilute each ligand 40-fold in **1x L3MBTL1 Assay Buffer**. Add 5 µL of diluted **Methylated MBT Ligand 1** to each well designated as "Positive Control" and "Test Inhibitor". Add 5 µL of diluted **Non-Methylated MBT Ligand 1** to the wells labeled as "Substrate Control".

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** info@bpsbioscience.com

- 7) Thaw **L3MBTL1** protein on ice. Upon first thaw, briefly spin tube containing protein to recover the full contents of the tube. Aliquot **L3MBTL1** protein into single-use aliquots. Store remaining undiluted **L3MBTL1** in aliquots at  $-80^{\circ}\text{C}$  immediately. *Note: L3MBTL1 is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 8) Dilute **L3MBTL1** in **1x L3MBTL1 Assay Buffer** to  $3.5\text{ ng}/\mu\text{l}$  ( $10.5\text{ ng}/\text{reaction}$ ). Initiate reaction by adding  $3\ \mu\text{l}$  of diluted **L3MBTL1** to wells designated for the "Substrate Control" "Positive Control", and "Test Inhibitor". Discard any remaining diluted L3MBTL1 protein after use.
- 9) Incubate at room temperature for 1 hour.
- 10) Read the fluorescent intensity in a microtiter-plate reader capable of TR-FRET.

#### Instrument Settings

Reading Mode	Time Resolved
Excitation Wavelength	$340\pm 20\text{ nm}$
Emission Wavelength	$620\pm 10\text{ nm}$
Lag Time	$60\ \mu\text{s}$
Integration Time	$500\ \mu\text{s}$
Excitation Wavelength	$340\pm 20\text{ nm}$
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).

When percentage activity is calculated, the FRET value from the substrate control can be set as zero percent activity and the FRET value from the positive control can be set as one hundred percent activity.

$$\% \text{ Activity} = \frac{\text{FRET}_s - \text{FRET}_{\text{Sub}}}{\text{FRET}_p - \text{FRET}_{\text{Sub}}} \times 100\%$$

Where  $\text{FRET}_s$  = Sample FRET,  $\text{FRET}_{\text{Sub}}$  = Substrate control FRET, and  $\text{FRET}_p$  = Positive control FRET.

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

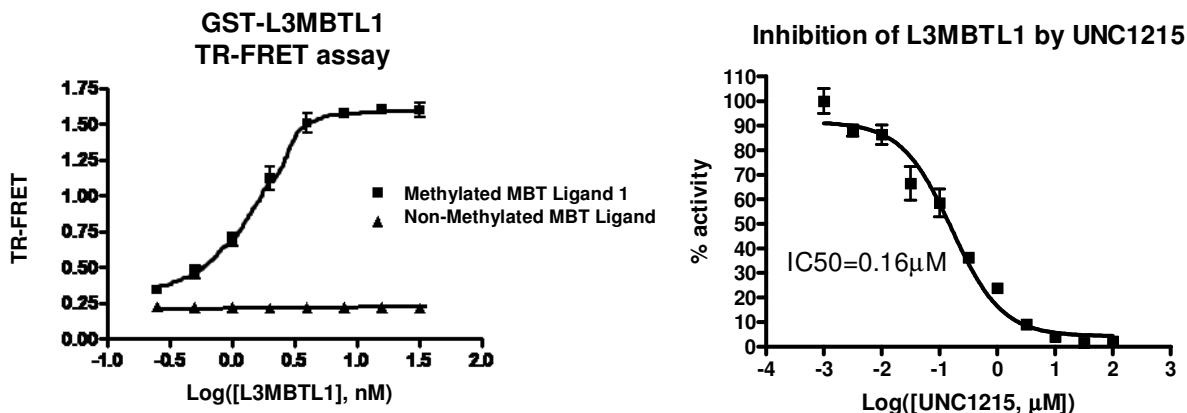
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
 San Diego, CA 92121  
 Tel: 1.858.829.3082  
 Fax: 1.858.481.8694  
 Email: info@bpsbioscience.com

**EXAMPLE OF ASSAY RESULTS:**



Interaction of L3MBTL1 (BPS Bioscience #55000) with peptide ligands, assayed using the L3MBTL1 TR-FRET Assay Kit (#55200). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com)*

**RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
L3MBTL1, GST-tag	55000	100 μg
L3MBTL1, His-tag	55002	100 μg
L3MBTL1 Inhibitor Screening Kit	55100	384 rxns.
L3MBTL2, His-tag	55018	100 μg
L3MBTL3, GST-tag	55013	100 μg
L3MBTL3, His-tag	55014	100 μg
UHRF1 (2-793), His-FLAG tag	55001	50 μg
UHRF1 (108-286), His-tag	55004	100 μg
UHRF1 (108-286), GST-tag	55003	100 μg
CBX1, GST-tag	55009	100 μg
CBX2, GST-tag	55011	100 μg
CBX7, His-tag	55017	100 μg
CHD2, GST-tag	55005	100 μg
CDY1, GST-tag	55007	100 μg
UNC1215	27404	1 mg
UNC1215	27405	5 mg
UNC1215	27406	10 mg

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)