



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
BCL2A1 TR-FRET Assay Kit
Catalog #79601
Size: 384 reactions

DESCRIPTION: The BCL2A1 TR-FRET Assay Kit is designed to measure the inhibition of BCL2A1 (BFL1) binding to its ligand 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, BCL2A1, peptide ligand, and an inhibitor is incubated for 2 hours. Then, the fluorescence intensity is measured using a fluorescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
100077	BCL2A1	20 µg	-80°C	(Avoid freeze/thaw cycles!)
	BCL2 Peptide Ligand 2	400 rxns	-80°C	
30017	Anti-His Tb-labeled donor	2 x 10 µl	-20°C	
	Dye-labeled acceptor	2 x 10 µl	-20°C	
	3x MCL TR-FRET Assay Buffer*	4 ml	-20°C	
	White, nonbinding, low volume, microtiter plate	1	Room temp.	

* Add 30 µl of 0.5M DTT before use.

MATERIALS REQUIRED BUT NOT SUPPLIED:

Dithiothreitol (DTT, 0.5 M)
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: Kvensakul, M., Hinds, MG. *Cell Death and Disease* 2013; **4**: e909.

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



6042 Cornerstone Court W, Ste 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 duplicate.

- 1) Add 30 μ l of 0.5M DTT to **3x MCL TR-FRET Assay Buffer** before use. Dilute one part **3x MCL TR-FRET Assay Buffer** with 2 parts distilled water (3-fold dilution) to make **1x MCL TR-FRET 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-labeled donor** and **Dye-labeled acceptor** 80-fold in **1x MCL TR-FRET Assay Buffer**. Make only sufficient quantities needed for the assay; store remaining stock solution in aliquots at -20°C .
- 3) Add 4 μ l of diluted **Anti-His Tb-labeled donor** and 4 μ l of diluted **Dye-labeled acceptor** to each well designated "Test Inhibitor," "Negative Control," and "Positive Control."
- 4) Add 4 μ l of inhibitor solution to each well designated "Test Inhibitor." Add 4 μ l of the same solution without inhibitor (inhibitor buffer) to the wells labeled "Negative Control" and "Positive Control."

	Negative Control*	Positive Control	Test Inhibitor
Anti-His Tb-labeled donor	4 μ l	4 μ l	4 μ l
Dye-labeled acceptor	4 μ l	4 μ l	4 μ l
Test Inhibitor	-	-	4 μ l
Inhibitor Buffer (no inhibitor)	4 μ l	4 μ l	-
1x MCL TR-FRET Assay Buffer	4 μ l	-	-
BCL2 Peptide Ligand 2	-	4 μ l	4 μ l
BCL2A1 (10 ng/ μ l)	4 μ l	4 μ l	4 μ l
Total	20 μl	20 μl	20 μl

- 5) Resuspend **BCL2 Peptide Ligand 2** in 1600 μ l of **1x MCL TR-FRET Assay Buffer**. Aliquot **BCL2 Peptide Ligand 2** into single-use aliquots. Store remaining ligand at -80°C immediately. *Note: Ligand is sensitive to freeze/thaw cycles. Do not re-use thawed aliquots.*
- 6) Add 4 μ l of resuspended **BCL2 Peptide Ligand 2** to each well designated as "Positive Control" and "Test Inhibitor." Add 4 μ l of **1x MCL TR-FRET Assay Buffer** to the wells labeled as "Negative Control."
- 7) Thaw **BCL2A1** protein on ice. Upon first thaw, briefly spin tube containing protein to recover the full contents of the tube. Aliquot **BCL2A1** protein into single-use aliquots. Store remaining undiluted aliquots at -80°C immediately. *Note: BCL2A1 is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*

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



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

- 8) Dilute **BCL2A1** in **1x MCL TR-FRET Assay Buffer** to 10 ng/μl (40 ng/reaction). Initiate reaction by adding 4 μl of diluted **BCL2A1** to wells designated for the “Negative Control,” “Positive Control,” and “Test Inhibitor.” Discard any remaining diluted **BCL2A1** protein after use.
- 9) Incubate at room temperature for 2 hours.
- 10) Read the fluorescent intensity in a microtiter-plate reader capable of TR-FRET.

Instrument Settings

Reading Mode	Time Resolved
Excitation Wavelength	340±20 nm
Emission Wavelength	620±10 nm
Lag Time	60 μs
Integration Time	500 μs
Excitation Wavelength	340±20 nm
Emission Wavelength	665±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).

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

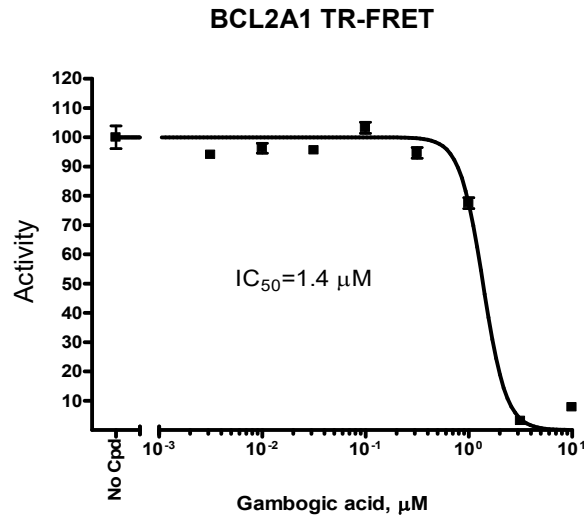
$$\% \text{ Activity} = \frac{FRET_S - FRET_{neg}}{FRET_P - FRET_{neg}} \times 100\%$$

Where $FRET_S$ = Sample FRET, $FRET_{Neg}$ = Negative control FRET, and $FRET_P$ = 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 **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com

EXAMPLE OF ASSAY RESULTS:



Inhibition of BCL2A1 by Gambogic Acid, measured using the *BCL2A1 TR-FRET Assay Kit*, BPS Bioscience #79601. 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>
BCL2A1	100077	100 μg
BCL2L2, His-Tag	100091	100 μg
BCL2L10, His-Tag	100080	100 μg
Bcl-2, His-tag	50272	100 μg
Bcl-xL, His-tag	50273	100 μg
MCL1, His-Tag	40742	100 μg
MCL-1 TR-FRET Assay Kit	79506	384 rxns.
BCL-2 TR-FRET Assay Kit	50222	384 rxns.
BCL-2L2 TR-FRET Assay Kit	79587	384 rxns.
BCL-xL TR-FRET Assay Kit	50223	384 rxns.
Caspase-3	80500	50 μg
Caspase-6	80113	50 μg
Caspase-7	70000	50 μg
Caspase-8	80114	50 μg
Caspase-9	80115	50 μg
Caspase-3 Homogeneous Assay Kit	80700	96 rxns.
Caspase-6 Homogeneous Assay Kit	80703	96 rxns.
Caspase-7 Homogeneous Assay Kit	80701	96 rxns.
Caspase-8 Homogeneous Assay Kit	80704	96 rxns.

Note: Anti-His Tb-labeled donor and dye-labeled acceptor are products of Cisbio Bioassays.

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