



6405 Mira Mesa Blvd. Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet
Cereblon Ubiquitination Homogeneous Assay Kit
Catalog #79881
Size: 384 reactions

BACKGROUND:

Cereblon (CRBN), a substrate receptor of Cullin-RING ligase 4 (CRL4), is a primary target of thalidomide and immunomodulatory imide drugs (IMiDs). CRBN binders are widely used in the design of heterobifunctional molecules called proteolysis-targeting chimeras (PROTACs) for degradation of the protein of interest. Several proteins, such as BRD4, CDK9, or Tau, can be successfully degraded by CRBN-based PROTACs. CRBN complex, CRL4^{CRBN}, is a unique E3 ubiquitin ligase because its substrate selectivity is altered by various ligands including IMiDs. IMiDs induce degradation of neosubstrates, such as Ikaros or CK1a. CRL4^{CRBN} also promotes auto-ubiquitination of CRBN, which can be significantly inhibited by thalidomide.

DESCRIPTION:

The *Cereblon Ubiquitination Homogeneous Assay Kit* is designed to measure Cereblon auto-ubiquitination activity in a homogeneous 384 reaction format. It utilizes biotin-labeled ubiquitin and proper A-screen beads to complete the pairing. This homogeneous assay requires no time-consuming washing steps, making it especially suitable for high throughput screening applications.

COMPONENTS:

Catalog #	Component	Amount	Storage	
100402	UBE1 (E1)	4 µg	-80°C	Avoid freeze/ thaw cycles!
80313	UBCH5C (E2)	40 µg	-80°C	
100405	Human Cereblon Complex (E3)	5 µg	-80°C	
	Biotin-Ubiquitin	400 µl	-80°C	
	ATP (40 mM)	20 µl	-80°C	
	U2 Assay Buffer	2 x 10 ml	-80°C	
	100 mM (-)-Thalidomide	15 µl	-80°C	

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



6405 Mira Mesa Blvd. Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

AlphaLISA anti-FLAG acceptor beads, 5 mg/ml (PerkinElmer #AL112C)
AlphaScreen Streptavidin-conjugated donor beads, 5 mg/ml (PerkinElmer #6760002S)
Optiplate 384 (PerkinElmer #6007290)
AlphaScreen microplate reader
Adjustable micropipettor and sterile tips

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

CONTRAINDICATIONS: Green and blue dyes that absorb light in the AlphaScreen signal emission range (520-620 nm), such as Trypan Blue. Avoid the use of the potent singlet oxygen quenchers such as sodium azide (NaN_3) or metal ions (Fe^{2+} , Fe^{3+} , Cu^{2+} , Zn^{2+} and Ni^{2+}). The presence of culture medium RPMI 1640 at >1% leads to signal reduction due to the presence of excess biotin and iron in this medium. MEM, which lacks these components, does not affect AlphaScreen assays.

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

REFERENCES:

- 1) Chamberlain PP, Hamann LG. *Nat Chem Biol.* 2019 Oct;**15**(10):937-944.
- 2) Chen, Y.-A., Peng, Y.-J., Hu, M.-C., Huang, J.-J., *et al.* *Scientific Reports* 2015; **5** (1).

SAFETY WARNING: Thalidomide is known to cause severe birth defects in humans. **It is very important to use all appropriate precautions when handling this compound.**

ASSAY PROTOCOL

All samples and controls should be tested at least in duplicate. All incubations at room temperature are performed with slow shaking on a rotator platform.

Step 1:

- 1) Thaw UBE1, UBCH5C, Cereblon, Biotin-Ubiquitin, ATP, and U2 Assay Buffer on ice. Aliquot each protein, assay buffer, and ATP into single-use aliquots and stored at -80°C immediately. *Note: UBE1, UBCH5C, Cereblon, Biotin-Ubiquitin, and assay buffer are sensitive to freeze/thaw cycles. Avoid multiple freeze-thaw cycles.*

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



6405 Mira Mesa Blvd. Suite 100
 San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 2) Carefully calculate the amount of proteins needed. Prepare appropriate amounts of diluted proteins and reagents as described below:

Dilute the Cereblon in U2 Assay Buffer at 2 ng/μl;
 Dilute the UBE1 in U2 Assay Buffer at 9 ng/μl;
 Dilute the UBCH5C in U2 Assay Buffer at 105 ng/μl;
 Dilute the ATP in U2 Assay Buffer at 2 mM (20-fold);
 Keep the diluted reagents on ice until use.

- 3) Add 4 μl of diluted Cereblon to each well designated “Positive Control” and “Test Sample”. For the “Blank”, add 4 μl of U2 Assay Buffer.
- 4) Add 2 μl of inhibitor solution to each well designated “Test Sample”. For the “Positive Control” and “Blank”, add 2 μl of the same solution without inhibitor (Inhibitor buffer).
- 5) Preincubate Cereblon complex with the inhibitor(s) at room temperature for 30 min with slow shaking.
- 6) During the preincubation, prepare a master mixture using the diluted reagents prepared in step 2): N wells × (1 μl Biotin-Ubiquitin + 1 μl UBE1 + 1 μl UBCH5C + 1 μl ATP). After 30 minutes, start the reaction by adding 4 μl of master mixture to each well.
- 7) Incubate the reaction at 30°C for two hours. Cover the plate with a plate sealer if necessary.

	Blank	Positive Control	Test Sample
Cereblon	–	4 μl	4 μl
U2 Assay Buffer	4 μl	–	–
Test Inhibitor/Activator	–	–	2 μl
Inhibitor buffer*	2 μl	2 μl	–
Biotin-Ubiquitin	1 μl	1 μl	1 μl
UBE1	1 μl	1 μl	1 μl
UBCH5C	1 μl	1 μl	1 μl
ATP (2 mM)	1 μl	1 μl	1 μl
Total	10 μl	10 μl	10 μl

*Inhibitor buffer typically represents U2 Assay buffer with proper concentration of DMSO.

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



6405 Mira Mesa Blvd. Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Note: Protect your samples from direct exposure to light for steps 2 and 3!

Step 2:

Thaw U2 Assay Buffer on ice. Dilute anti-FLAG Acceptor beads (PerkinElmer #AL112C) 250-fold with U2 Assay Buffer. Add 10 μ l per well. Shake on a rotator platform for 30 minutes at room temperature.

Step 3:

1) Dilute Streptavidin-conjugated donor beads (PerkinElmer #6760002S) 125-fold with U2 Assay Buffer. Add 10 μ l per well. Shake on a rotator platform for 15-30 minutes at room temperature.

2) Read Alpha-counts. "Blank" value should be subtracted from all readings.

Example of Assay Results:

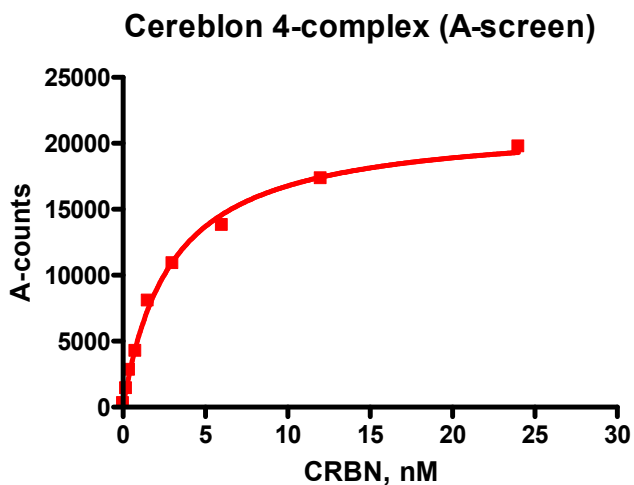


Figure 1: Titration of Cereblon ubiquitination using the *Cereblon Ubiquitination Homogeneous Assay Kit*, BPS Bioscience #79881. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com.

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

CEREBLON ubiquitination (A-screen)

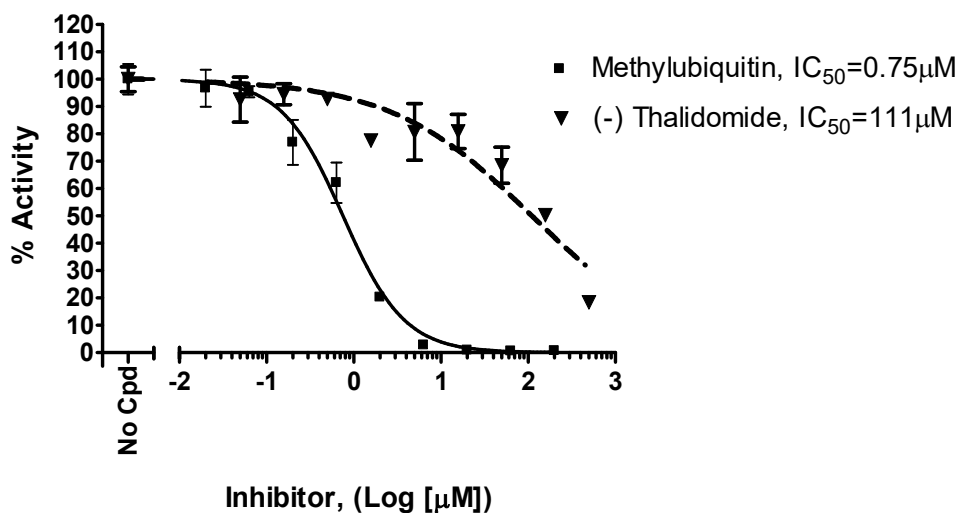


Figure 2: Inhibition of Cereblon ubiquitination by Methylated Ubiquitin or (-)Thalidomide, measured using the *Cereblon Ubiquitination Homogeneous Assay Kit*, BPS Bioscience #79881. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com.

RELATED PRODUCTS

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
CBL-B, GST-Tag (Human)	#80415	100 µg
CBL-B, His-Avi-Tag	#80414	100 µg
CBL-B, Biotin-labeled (Human)	#80412	50 µg
CBL-B (Y363F), Biotin-labeled (Human)	#80413	50 µg
UBE1 (UBA1), FLAG-tag	#80301	100 µg
UBCH5b	#80314	100 µg
Cereblon/DDB1/Cul4A/Rbx1 Complex	#100329-1	10 µg
PROTAC Optimization Kit for BET Bromodomain-Cereblon Binding	#79770	384 rxns.

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