

Data Sheet

BRD2 (339-459), GST-tag

Human, recombinant, N-terminal GST-tag
Catalog #: 31023
Lot #: 150311-1 **Conc.:** 3.2 mg/ml

Formulated in: 40 mM Tris-HCL, pH 8.0,
 110 mM NaCl, 2.2 mM KCl and 20%
 glycerol

Stability: At least 6 months at -80°C .
 Avoid freeze/thaw cycles. Storing diluted
 protein is not recommended, if necessary,
 use carrier protein (BSA 0.1 – 0.5%).

References:

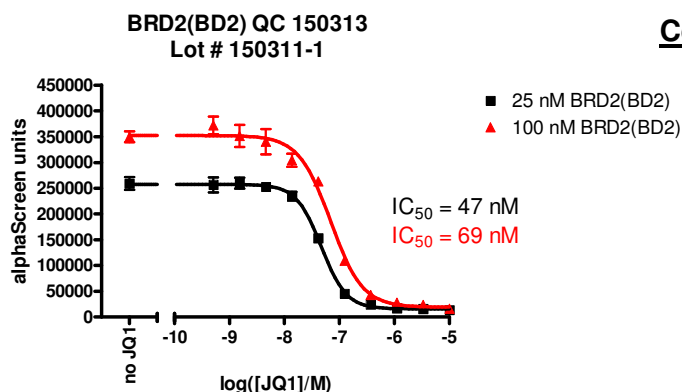
1. Umehara, T., *et al.*, *FEBS Lett.* 2010
 Sep 24; **584(18)**: 3901-3908.
2. Shang, E., *et al.*, *J Cell Biochem.* 2011
 Oct; **112(10)**: 2784-2793.

Description: Human Bromodomain
 containing 2, also known as BRD2, RING3
 or RNF3, amino-acids 339 – 459
 (GenBank Accession No. NM_005104)
 with N-terminal GST tag, MW = 41.0 kDa,
 expressed in an *E. coli* expression system.

Assay Conditions: Assay was performed
 using BRD2 BD2 Inhibitor Screening
 Assay Kit, Cat. #32522. BRD2 protein is
 incubated with biotinylated acetyl-histone
 peptide substrate (Cat. #33000) in BRD
 assay buffer (Cat. #33001) in a 10 μl
 reaction for 1 hour at RT. AlphaLISA®
 GSH acceptor beads (Perkin Elmer) are
 added, followed by AlphaScreen®
 Streptavidin-conjugated donor beads
 (Perkin Elmer). After 1 hour incubation at
 RT, Alphacounts are measured.

Application: Useful for the study of
 bromodomain binding assays, screening
 inhibitors, and selectivity profiling.

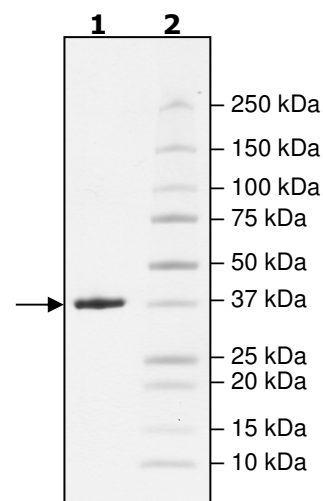
Quality Assurance



4-20% SDS-PAGE **Coomassie staining**

Lane 1:
 1.8 μg BRD2
Lane 2:
 Protein Marker

MW: 41 kDa
Purity: $\geq 95\%$



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

Please visit our website at: www.bpsbioscience.com