

Data Sheet

BRD2 (339-459)

Human, recombinant, N-terminal HIS tag

Catalog #: 31020

Lot #: 120207G Conc.: 0.9 mg/ml

Formulated in: 50 mM Tris-HCl, pH 8.0, 138 mM NaCl, 2.7 mM KCl, 0.05% Tween-20, and 10% glycerol.

Stability: >6 months at -80 °C

References:

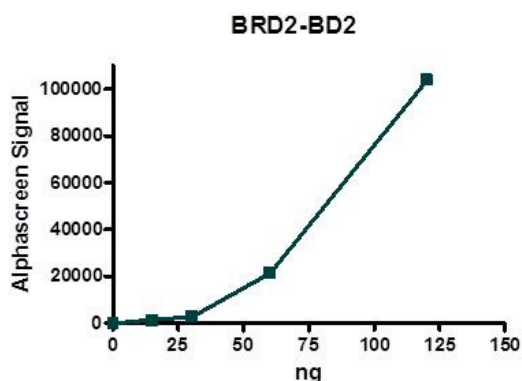
1. Umehara, T., *et al.*, *FEBS Lett.* 2010 Sep 24; **584(18)**: 3901-3908.
2. Shang, E., *et al.*, *J Cell Biochem.* 2011 Oct; **112(10)**: 2784-2793.

Description: Human Bromodomain containing 2, also known as BRD2, RING3 or RNF3, amino-acids 339 - 459 (GenBank Accession No. NM_005104) with N-terminal HIS tag, MW = 15.1 kDa, expressed in an *E. coli* expression system.

Assay Conditions: BRD2 is incubated with biotinylated acetyl-histone peptide substrate (BPS #33000) in BRD assay buffer (BPS #33001) in a 10 µl reaction for 1 hour at RT. AlphaLISA® Nickel Chelate 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

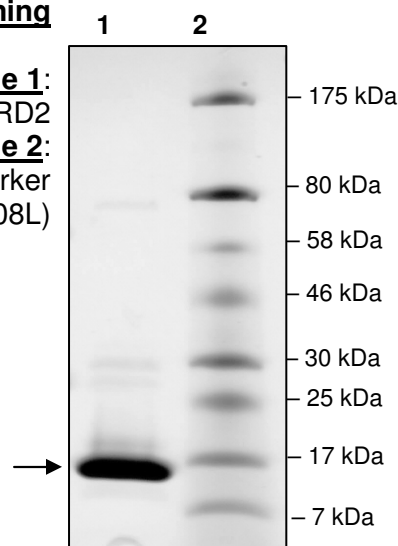


4-20% SDS-PAGE
Coomassie staining

Lane 1:
4.2 µg BRD2

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 15.1 kDa
Purity: ≥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