

Data Sheet

BRD2, BD1 and BD2 (65-459)

Human, recombinant, N-terminal His-tag

Catalog#: 31025

Lot#: 121130 **Conc.:** 1.42 mg/ml

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

Stability: >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 65 – 459, GenBank
Accession No. NM_005104, with N-
terminal His-tag, MW = 45 kDa,
expressed in an *E. coli* expression
system.

Application:

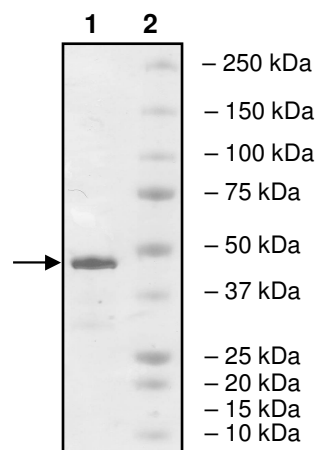
Useful for the study of bromodomain
binding assays, screening inhibitors, and
selectivity profiling.

Quality Assurance

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

Lane 1:
2 μg BRD2
Lane 2:
Protein Marker

MW: 45 kDa
Purity: $\geq 90\%$



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



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

Protein Sequence

MHHHHHEVSNPKKPGRVTNQLQYLHKVVMKALWKHQFAWPFRQPVDAVKLGLPDYH
KIIKQPMDMGTIKRRELENNYYWAASECMQDFNTMFTNCYIYNKPTDDIVLMAQTLEKIFLQ
KVASMPQEEQELVVTIPKNSHKKGAKLAALQGSVTSAHQVPAVSSVSHTALYTPPEIPT
TVLNIPHSVISSPLLKSLHSAGPPLLAVTAAPPAQPLAKKKGVKRKADTTTPTPTAILAPG
SPASPPGSLEPKAARLPPMRRESGRPIKPPRKDLPDSQQQHQSCKKGLSEQLKHCNGI
LKELLSKKHAAYAWPFYKPVDASALGLHDYHDIKHPMDLSTVKRKMENRDYRDAQEFA
ADVRLMFSNCYKYNPPDHDVVAMARKLQDVFEFRYAKMPDEPLE

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