

Data Sheet

UHRF1 (108-286)

Human, recombinant, N-terminal His tag

Catalog #: 55004

Lot #: 140205 Conc.: 0.6 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 10 mM β -mercaptoethanol, 250 mM imidazole, and 20% glycerol.

Stability: >6 months at -80°C

References:

1. Kofunato, Y., *et al.*, *Oncol Rep.* 2012 Dec; **28(6)**:1997-2002.
2. Jazirehi, A.R., *et al.*, *Epigenomics.* 2012 Jun; **4(3)**:251-252.

Description:

Tandem Tudor Domain region (a.a. 108-286) of human ubiquitin-like with PHD and ring finger domains 1 (UHRF1), GenBank Accession #NM_001048201, with N-terminal His tag, MW = 21.6 kDa, expressed in an *E. coli* cell expression system.

Application:

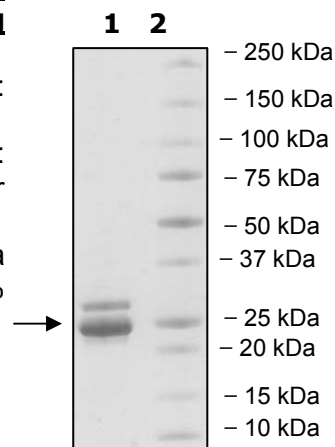
Useful for protein binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg UHRF1
Lane 2:
Protein Marker

MW: 21.6 kDa
Purity: $\geq 80\%$



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

Amino Acid Sequence:

MHHHHHHSTHGAAAETDSRPADEDMWDETELGLYKVNEYVDARDTNMGAWFEAQ
VVRVTRKAPSRDEPCSSSRPALEEDVIYHVKYDDYPENGVVQMNSRDVRRARARTIIK
WQDLEVGQVVMLNYPDNPKERGFWDYDAEISRKRETRTARELYANVVLGDDSLNDC
RIIFVDEVFKIERPGEG

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