

Data Sheet

A20 (TNFAIP3)

Human, recombinant, N-terminal His-tag

Catalog #: 80393

Lot #: 120924 Conc.: 3.8 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 250 mM Imidazole, and 20% Glycerol.

Stability: >6 months at -80 °C

References:

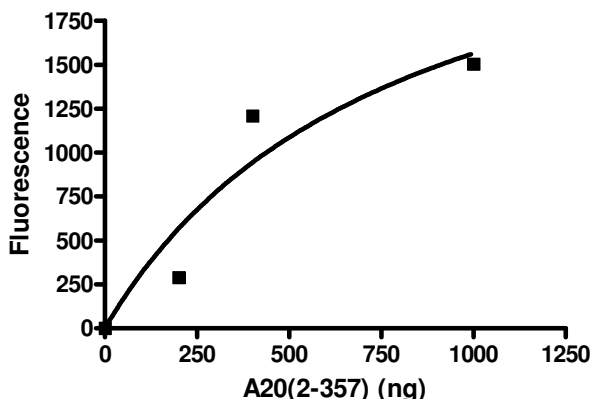
1. Shi CS, Kehrl JH. (2010). *Autophagy* **6(7)**:986-7.
2. Tavares, R.M., et al. (2010). *Immunity* **33(2)**:181-9.
3. Hymowitz, S.G., Wertz, I.E. (2010). *Nat. Rev. Cancer* **10(5)**:332-41.

Description: Human A20, also known as tumor necrosis factor- α induced protein 3 (TNFAIP3) or OTU domain-containing protein 7C (OTUD7C), GenBank Accession No. NM_006290, amino-acids 2-357 with an N-terminal His-tag, MW=42.8 kDa, expressed in an *E. coli* cell expression system.

Specific Activity: 10.5 pmol/min/ μ g
 Assay conditions: Reaction was performed in 50 mM Tris, pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and A20. Reaction was incubated at 37°C for 15 min and fluorescent signal was measured at excitation = 340 nm, and emission = 460 nm.

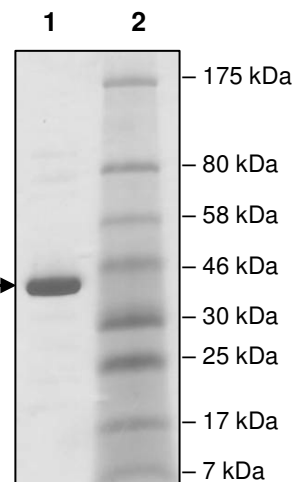
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 3 μ g A20
Lane 2:
 Protein Marker
MW: 42.8 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



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

Sequence:

MHHHHHHAEQVLPQALYLSNMRKAVKIRERTPEDIFKPTNGIIHHFKTMHRYTLEMFR
TCQFCPQFREIIHKALIDRNIQATLESQKLNWCREVRKLVALKTNGDGNCLMHATSQ
YMWGVQDSDLVLRKALFSTLKETDTRNFKFRWQLESLSQEFVETGLCYDTRNWND
EWDNLIKMASTDTPMARSGLQYNSLEEIHIFVLCNILRRPIIVISDKMLRSLESGSNFAP
LKVGGIYLPLHWPQAQECYRYPIVLGYDSHHFVPLVTLKDSGPEIRAVPLVNRDRGRFE
DLKVHFLTPENEMKEKLLKEYLMVIEIPVQGWDHGTTHLINAACLDEANLPKEINLVD
DYFELVQHEYKKWQ

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