

Data Sheet

Ataxin-3

Human, recombinant, N-terminal His-tag

Catalog #: 80399

Lot #:101019

Conc.: 1.18 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM imidazole, 10% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

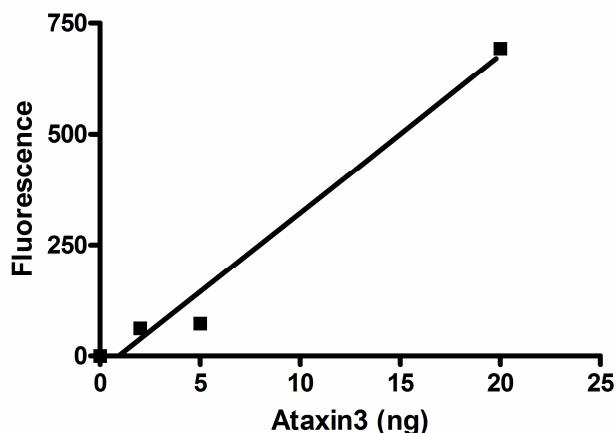
1. do Carmo Costa, M., *et al.* (2010). *PLoS One* **5(7)**:e11728.
2. Jung, J., *et al.* (2009). *Hum Mol Genet.* **18(24)**:4843-52.
3. Macedo-Ribeiro, S., *et al.* (2009). *PLoS One* **4(6)**: e5834.

Description: Human Ataxin-3 (ATX3 or ATXN3), also known as Josephin (JOS), Machado-Joseph disease protein 1 (MJD1) or Spinocerebellar ataxia type 3 (SCA3), GenBank Accession No. NM_004993, amino-acids 2-362(end) with an N-terminal His-tag, MW=42.1 kDa, expressed in *E. coli*.

Activity Assay: Reaction was performed in 50 mM Tris, pH 7.4, 1 mM DTT, 0.5 mM EDTA, 500 nM Ub-AMC, and Ataxin-3 (0-0.4 ng/μL) in a 50 μL reaction. Reaction was incubated at 37°C for 15 min and fluorescent signal was measured at excitation = 340 nm, emission = 460 nm.

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

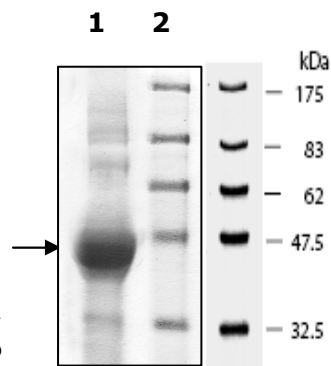
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 10 μg Ataxin-3
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 42 kDa.
Purity: ≥60%



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