

Data Sheet

Ubiquitin, His-Avi-Tag, Biotin Labeled

Human, Recombinant, N-terminal His-Avi-Tag

Catalog #: 11236

Lot #: 181207-2

Conc: 0.70 mg/ml

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

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

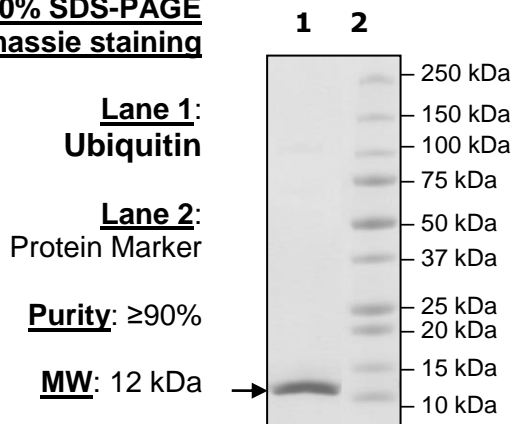
1. Bodine, S., *et al.*, *Science*. 2001; **294(5547)**: 1704-1708.
2. Bence, N., *et al.*, *Science*. 2001; **292(5521)**: 1552-1555.

Description:

Human Ubiquitin, also known as UBB, UBC, UBA52, and RPS27A. GenBank Accession No. NM_021009, a.a. 2-76, with N-terminal His-Avi-Tag™, biotin labeled, expressed in an Sf9 expression system. This protein is enzymatically biotinylated using Avi-Tag technology. MW = 12 kDa.

Quality Assurance

4-20% SDS-PAGE Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.481.8694**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com