

Data Sheet

USP9X, His-tag, FLAG-tag

Human, recombinant, N-terminal His-tag,
C-terminal FLAG-tag

Catalog #: 100188

Lot #: 190409

Conc.: 0.63 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, 20% Glycerol and 100 µg/mL FLAG
peptide.

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

References:

1. Schwickart, M., *et al.*, *Nature*. 2010,
Jan; **463**:103-107.
2. Dupont, S., *et al.*, *Cell*. 2009, Jan
1(9):13-14.

Description: Human USP9X also known
as Ubiquitin Specific Peptidase 9 X-
Linked, FAM, DFFRX, Ubiquitin
Thioesterase FAF-X, MRXS99 and
Deubiquitinating Enzyme FAF-X,
GenBank Accession No.
NM_001039591, a.a. 1559-1995 with N-
terminal His-tag and C-terminal FLAG-
tag, expressed HEK293 expression
system. MW = 53 kDa.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

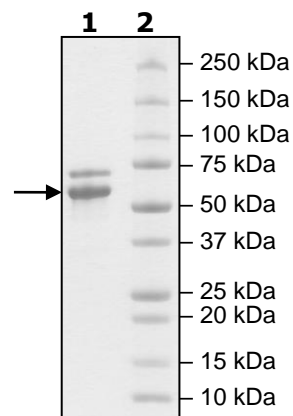
2 µg USP9X

Lane 2:

Protein Marker

MW: 53 kDa

Purity: ≥72%



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** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com