

Data Sheet

FKBP12, His-tag

Human, recombinant, N-terminal His-tag

Catalog #: 40550

Lot#: 90812 **Conc.:** 3.5 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM EGTA, 0.1 mM EDTA, 0.1 mM PMSF, 50% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Knappe, T.A. et al., J. Mol. Biol. 368 (5), 1458-1468 (2007).
2. Wear, M.A. et al., Protein Expr. Purif. 51 (1), 80-95 (2007).

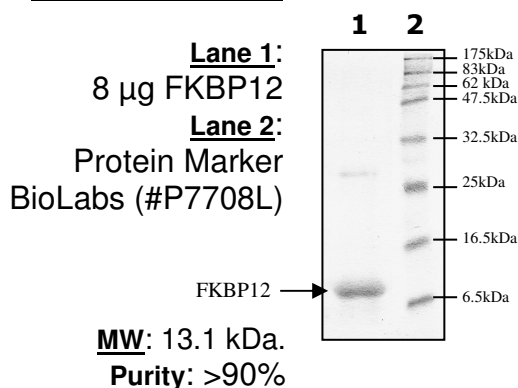
Description: Human FKBP12 (FK506 binding protein 1A, transcript variant 12A) (GenBank Accession No. NM_054014), full length with N-terminal His tag, MW= 13.1 kDa, expressed in E. Coli expression system.

Application:

Useful for testing rapamycin inhibition of mTOR.

Quality Assurance

**14% 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 or **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