

Data Sheet

TUT4, FLAG-tag, Biotin-Labeled

Human, Recombinant, C-terminal FLAG-Avi-tag
Catalog #: 71258

Lot #: 160819-D **Conc.:** 0.87 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20 and 20% 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%).

Description: Human TUT4, also known as
ZCCHC11, GenBank Accession number
NM_001009881, a.a. 212-1313 with C-
terminal FLAG-Avi-Tag. MW = 127 kDa,
expressed in an HEK293 expression
system. This protein is enzymatically
biotinylated using Avi-Tag™ technology.
Biotinylation is confirmed to be $\geq 90\%$

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

References:

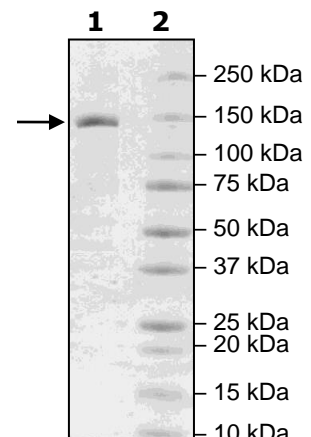
1. Hagan, J. *et al.*, *Nature Structural & Molecular Biology*. 2009 Aug **16**: 1021-1025
2. Jones, M. *et al.*, *Nature Cell Biology*. 2009 Aug; **11**:1157-1163

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg TUT4
Lane 2:
Protein Marker

MW: 127 kDa
Purity: $\geq 89\%$



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