

Data Sheet

ADAR1, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 100472

Lot #: 191206-1

Conc.: 0.24 mg/mL

Formulated in: 50 mM Tris-HCl, pH 8.0,
750 mM NaCl, 0.01% Triton X-100, 10%
glycerol and 100 µg/ml FLAG peptide.

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

References:

1. Wang, Q., *et al.*, *Science*. 2000;
290(5497):1765-1768.
2. Wong, S.K., *et al.*, *RNA*. 2001; **7:**846-
858

Description: Human adenosine deaminase, RNA-specific 1 (ADAR1) protein, also known as DSH, and DRADA, GenBank Accession No. NM_001111, a.a 2-1227(end) with an N-terminal FLAG-tag, expressed in a HEK293 cell expression system. MW =137 kDa.

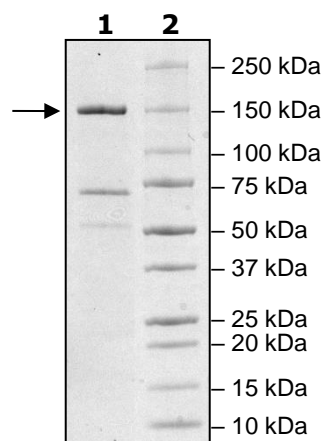
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2µg ADAR1
Lane 2:
Protein Marker

MW: 137 kDa

Purity: ≥70%



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