

Data Sheet

RbAp48

Human, recombinant, N-terminal His-tag

Catalog #: 50297

Lot#: 111011 **Conc.:** 0.48 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 200 mM
Imidazole, 20% Glycerol, and 3mM DTT.

Stability: >6 months at -80°C

References:

1. Varambally, S. *et al. Science* **322 (5908)**, 1695-1699 (2008).
2. Rakotobe, D. *et al., Virol. J.* **5**, 32 (2008).

Description:

Human RbAp48 (GenBank Accession No. NM_005610), a.a. 2-end, with N-terminal His-tag, MW = 48 kDa, expressed in a baculovirus infected Sf9 cell expression system.

Application:

Useful for western blot analysis.

Quality Assurance

10% SDS-PAGE
Coomassie staining

Lane 1:

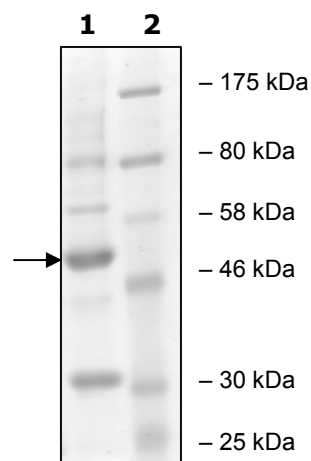
1.7 µg RbAp48

Lane 2:

Protein Marker
BioLabs (#P7708L)

MW: 48 kDa

Purity: ≥ 46%



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