

Data Sheet

AXL, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 100174

Lot#: 180907 **Conc.:** 0.77 mg/ml

Description: Human AXL, also known as Tyrosine-protein kinase receptor UFO, and JTK11, GenBank Accession No. NM_021913, a.a. 473-894 (end) with N-terminal FLAG-Tag, MW= 49 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 100ug/ml flag

Stability: At least 6 months at -80°C

References:

1. Sawabu T. et al., Mol Carcinog. 2007 Feb;46(2):155-64.
2. Vajkoczy P. et al., Proc Natl Acad Sci U S A. 2006 Apr 11;103(15):5799-804.

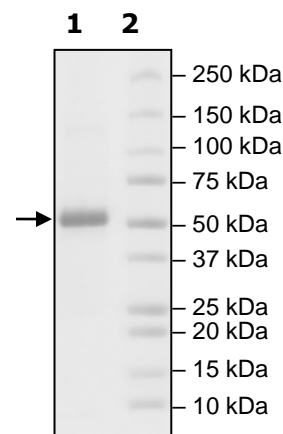
Quality Assurance

4-20% SDS-PAGE Coomassie Staining

Lane 1:
AXL

Lane 2:
Protein Marker

Purity: >90%
MW: 49 kDa



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 **1.858.259.5734**

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

Please visit our website at: www.bpsbioscience.com