

Catalogue #	Aliquot Size
A34-12BG -05	5 µg
A34-12BG -10	10 µg

# AXL (R499C), Active

Recombinant protein expressed in Sf9 cells

#### Catalog # A34-12BG Lot # J570-3

### **Product Description**

Recombinant human AXL (R499C) (473-end) was expressed by baculovirus in Sf9 insect cells using an Nterminal GST tag. The Axl gene accession number is <u>NM 021913</u>.

### **Gene Aliases**

UFO, JTK11

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### **Storage and Stability**

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

## **Scientific Background**

AXL(R499C) is an arginine to cysteine mutation at 499 aa in AXL and appears as a somatic mutation in gastric adenocarcinoma samples. AXL is a transforming gene isolated from primary human myeloid leukemia cells, encodes a novel receptor tyrosine kinase which is a member of the TAM (TYRO3/AXL/MER) receptor tyrosine kinase subfamily (1). AXL represents a unique structure of the extracellular region that juxtaposes IgL and FNIII repeats and AXL transduces signals from the extracellular matrix into the cytoplasm by binding growth factors like vitamin K-dependent protein growth-arrest-specific gene 6 (2).

### References

- 1. O'Bryan, J. P.et.al: Axl, a transforming gene isolated from primary human myeloid leukemia cells, encodes a novel receptor tyrosine kinase. Molec. Cell. Biol. 11: 5016-5031, 1991.
- 2. Varnum, B. C.et.al: Axl receptor tyrosine kinase stimulated by the vitamin K-dependent protein encoded by growtharrest-specific gene 6. Nature 373: 623-626, 1995.

## **Specific Activity**



The specific activity of AXL (R499C) was determined to be **3 nmol** /min/mg as per activity assay protocol.

## Purity



The purity of AXL (R499C) was determined to be >85% by densitometry, approx. MW 76 kDa.

# AXL (R499C), Active

Recombinant human protein expressed in Sf9 cells

Catalog Number Specific Activity Specific Lot Number Purity Concentration

Storage & Shipping

A34-12BG 3 nmol/min/mg J570-3 >85% 0.1 µg/µl 1yr at -70°C from date of shipment Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: <u>orders@signalchem.com</u> <u>www.signalchem.com</u>

# Activity Assay Protocol

**Reaction Components** 

### Active Kinase (Catalog #: A34-12BG)

Active AXL (R499C)  $(0.1\mu g/\mu I)$  diluted with Kinase Dilution Buffer III (Catalog #: K23-09) and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active AXL (R499C) for optimal results).

### Kinase Dilution Buffer III (Catalog #: K23-09)

Kinase Assay Buffer I (Catalog #: K01-09) diluted at a 1:4 ratio (5X dilution) with 50ng/µl BSA solution.

### Kinase Assay Buffer I (Catalog #: K01-09)

Buffer components: 25mM MOPS, pH 7. 2, 12.5mM  $\beta$ -glycerol-phosphate, 25mM MgC1<sub>2</sub>, 5mM EGTA, 2mM EDTA. Add 0.25mM DTT to Kinase Assay Buffer prior to use.

### [<sup>33</sup>P]-ATP Assay Cocktail

Prepare 250µM [<sup>33</sup>P]-ATP Assay Cocktail in a designated radioactive working area by adding the following components: 150µl of 10mM ATP Stock Solution (Catalog #: **A50-09**), 100µl [<sup>33P</sup>]-ATP (1mCi/100µl), 5.75ml of Kinase Assay Buffer I (Catalog #: **K01-09**). Store 1ml aliquots at -20°C.

### 10mM ATP Stock Solution (Catalog #: A50-09)

Prepare ATP stock solution by dissolving 55mg of ATP in 10ml of Kinase Assay Buffer I (Catalog #: K01-09). Store 200 $\mu$ l aliquots at -20°C.

Substrate (Catalog #: A16-58)

AXLtide peptide substrate (KKSRGDYMTMQIG) diluted in distilled  $H_2O$  to a final concentration of 1mg/ml.

#### Assay Protocol

- Step 1. Thaw [<sup>33</sup>P]-ATP Assay Cocktail in shielded container in a designated radioactive working area.
- Step 2. Thaw the Active AXL (R499C), Kinase Assay Buffer, Substrate and Kinase Dilution Buffer on ice.
- **Step 3.** In a pre-cooled microfuge tube, add the following reaction components bringing the initial reaction volume up to 20μl:
  - Component 1. 10µl of diluted Active AXL(R499C) (Catalog # A34-12BG)
  - Component 2. 5µl of 1mg/ml stock solution of substrate (Catalog # A16-58)
  - Component 3. 5µl distilled H<sub>2</sub>O (4°C)
- Step 4. Set up the blank control as outlined in step 3, excluding the addition of the substrate. Replace the substrate with an equal volume of distilled H<sub>2</sub>O.
- Step 5. Initiate the reaction by the addition of 5 μl [<sup>33</sup>P]-ATP Assay Cocktail bringing the final volume up to 25μl and incubate the mixture in a water bath at 30°C for 15 minutes.
- **Step 6.** After the 15 minute incubation period, terminate the reaction by spotting 20 μl of the reaction mixture onto individual pre-cut strips of phosphocellulose P81 paper.
- Step 7. Air dry the pre-cut P81 strip and sequentially wash in a 1% phosphoric acid solution (dilute 10ml of phosphoric acid and make a 1L solution with distilled H<sub>2</sub>O) with constant gentle stirring. It is recommended that the strips be washed a total of 3 intervals for approximately 10 minutes each.
- Step 8. Count the radioactivity on the P81 paper in the presence of scintillation fluid in a scintillation counter.
- Step 9. Determine the corrected cpm by removing the blank control value (see Step 4) for each sample and calculate the kinase specific activity as outlined below.

## Calculation of [P<sup>33</sup>]-ATP Specific Activity (SA) (cpm/pmol)

Specific activity (SA) = cpm for 5 µl [<sup>33</sup>P]-ATP / pmoles of ATP (in 5 µl of a 250 µM ATP stock solution, i.e., 1250 pmoles)

### Kinase Specific Activity (SA) (pmol/min/µg or nmol/min/mg)

Corrected cpm from reaction / [(SA of <sup>33</sup>P-ATP in cpm/pmol)\*(Reaction time in min)\*(Enzyme amount in µg or mg)]\*[(Reaction Volume) / (Spot Volume)]

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: <u>orders@signalchem.com</u> www.signalchem.com

#### FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.