

PRODUCT INFORMATION



Axl Extracellular Domain (human, recombinant)

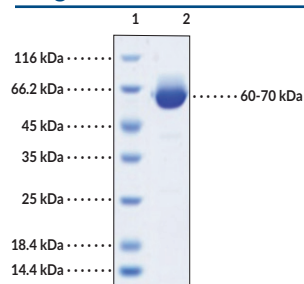
Item No. 44185

Overview and Properties

Synonym: UFO
Source: Active recombinant human C-terminal His-tagged Axl extracellular domain expressed in HEK293 cells
Amino Acids: 33-449
Uniprot No.: P30530
Molecular Weight: 46.5 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥98% estimated by SDS-PAGE
Supplied in: Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay

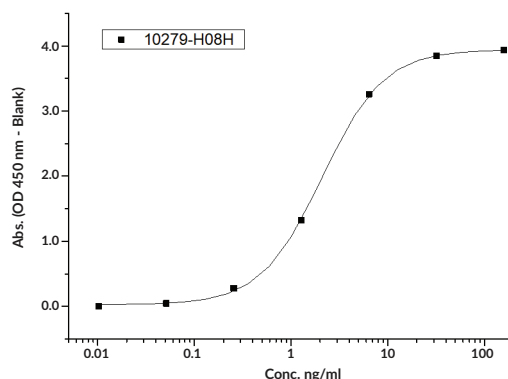
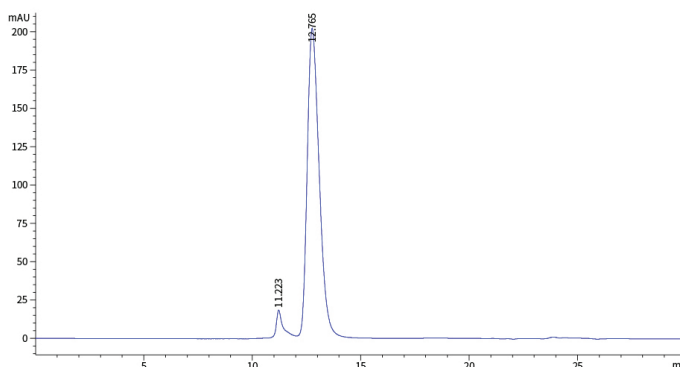
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images

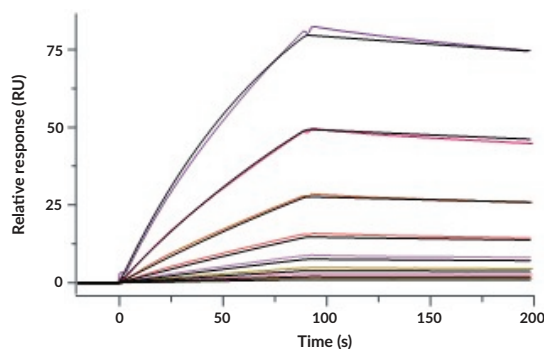


Lane 1: MW Markers
Lane 2: Axl Extracellular Domain

SDS-PAGE Analysis of ACTRIIB Extracellular Domain. This protein has a calculated molecular weight of 46.5 kDa. It has an apparent molecular weight of approximately 60-70 kDa by SDS-PAGE under reducing conditions due to glycosylation.



Label Recombinant Mouse GAS6 Protein (His Tag) with biotin. Immobilized Axl Extracellular Domain (human, recombinant) at 0.25 μg/ml (100 μl/well) can bind biotinylated Recombinant Mouse GAS6 Protein, the EC₅₀ is 1.3-3.9 ng/ml (routinely tested).



Axl Extracellular Domain (human, recombinant) (Item No. 44185) immobilized on CM5 chip, can bind recombinant human GAS6 protein (His-tag) with an affinity constant of 5.700 nM as determined in an SPR assay (Biacore 8K) (QC tested).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/20/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Axl is a receptor tyrosine kinase and member of the TAM kinase family.^{1,2} It is composed of two immunoglobulin-like (Ig-like) domains, two fibronectin-like III domains, an extracellular cleavage site, a transmembrane region, and a C-terminal kinase domain.¹ It is ubiquitously expressed and localizes to the cell membrane.^{3,4} After homodimerization, Axl is activated by growth arrest-specific protein 6 (GAS6) and is involved in promoting proliferation and hematopoietic lineage differentiation and negatively regulating antigen-presenting cell activation.^{1,2,5} The extracellular domain of Axl can be cleaved by metalloproteinases, releasing soluble Axl to bind and activate membrane-associated Axl.^{3,4} In cancer, Axl promotes metastasis, epithelial-to-mesenchymal transition (EMT), and survival.⁵ Increased expression of AXL is associated with shorter progression-free survival in patients with non-small cell lung cancer (NSCLC).⁶ Cayman's Axl Extracellular Domain (human, recombinant) protein can be used for binding assays. This protein consists of 428 amino acids, has a calculated molecular weight of 46.5 kDa, and a predicted N-terminus of Glu33 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 60-70 kDa due to glycosylation.

References

1. Graham, D.K., DeRyckere, D., Davies, K.D., *et al.* The TAM family: Phosphatidylserine sensing receptor tyrosine kinases gone awry in cancer. *Nat. Rev. Cancer* **14**(12), 769-785 (2014).
2. Axelrod, H. and Pienta, K.J. Axl as a mediator of cellular growth and survival. *Oncotarget*. **5**(19), 8818-8852 (2014).
3. Verma, A., Warner, S.L., Vankayalapati, H., *et al.* Targeting Axl and Mer kinases in cancer. *Mol. Cancer Ther.* **10**(10), 1763-1773 (2016).
4. Tondo, G., Perani, D., and Comi, C. TAM receptor pathways at the crossroads of neuroinflammation and neurodegeneration. *Dis. Markers* 2387614 (2019).
5. Myers, K.V., Amend, S.R., and Pienta, K.J. Targeting Tyro3, Axl and MerTK (TAM receptors): Implications for macrophages in the tumor microenvironment. *Mol. Cancer* **18**(1), 94 (2019).
6. Yoshimura, A., Yamada, T., Serizawa, M., *et al.* High levels of AXL expression in untreated EGFR-mutated non-small cell lung cancer negatively impacts the use of osimertinib. *Cancer Sci.* **4**(2), 606-618 (2023).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM