

# PRODUCT INFORMATION



## ALK2 Extracellular Domain (human, recombinant)

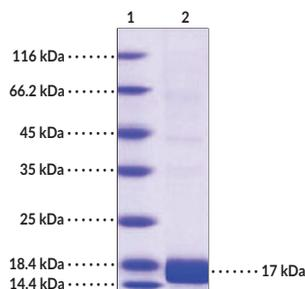
Item No. 44183

### Overview and Properties

<b>Synonyms:</b>	Activin Receptor-like Kinase 2, Activin Receptor Type I, ACTR-I, ACVR1, Serine/threonine-protein Kinase Receptor R1, TGF- $\beta$ Superfamily Receptor Type I, TSR-I
<b>Source:</b>	Recombinant human C-terminal His-tagged ALK2 extracellular domain expressed in insect cells
<b>Amino Acids:</b>	21-124
<b>Uniprot No.:</b>	Q04771
<b>Molecular Weight:</b>	12.8 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	$\geq 1$ year
<b>Purity:</b>	$\geq 93\%$ estimated by SDS-PAGE
<b>Supplied in:</b>	Sterile 20 mM Tris, pH 7.4, with 500 mM sodium chloride and 10% glycerol
<b>Endotoxin Testing:</b>	<1.0 EU/ $\mu$ g, determined by the LAL endotoxin assay
<b>Bioactivity:</b>	See figures for details

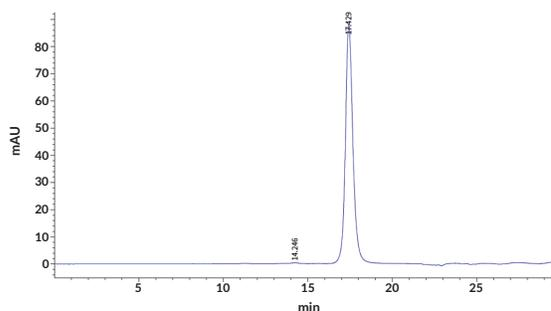
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: ALK2 Extracellular Domain

**SDS-PAGE Analysis of ALK2 Extracellular Domain.** This protein has a calculated molecular weight of 12.8 kDa. It has an apparent molecular weight of approximately 17 kDa by SDS-PAGE under reducing conditions due to glycosylation.



Size Exclusion Chromatography of purified ALK2 Extracellular Domain.

**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, 01/07/2026

**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

---

Activin receptor-like kinase 2 (ALK2) is a type I receptor kinase and member of the bone morphogenetic protein (BMP) receptor family.<sup>1,2</sup> It is composed of a ligand-binding extracellular domain, a transmembrane domain, a membrane-associated glycine/serine-rich (GS) domain, and an intracellular kinase domain. ALK2 is expressed primarily in the heart, liver, intestine, and kidneys but is found at low levels in the brain and lung.<sup>3</sup> Upon ligand binding, ALK2 is activated by type II BMP receptors via phosphorylation of the GS domain, which then activates SMAD or MAPK signaling.<sup>1,2</sup> SNPs in ACVR1, the gene encoding ALK2, are associated with fibrodysplasia ossificans progressiva (FOP), congenital heart defects, and diffuse intrinsic pontine glioma.<sup>4-6</sup> Cayman's ALK2 Extracellular Domain (human, recombinant) protein consists of 114 amino acids, has a calculated molecular weight of 12.8 kDa, and a predicted N-terminus of Met21 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 17 kDa due to glycosylation.

## References

---

1. Shi, F., Gao, J., Zou, J., *et al.* Targeting heterotopic ossification by inhibiting activin receptor-like kinase 2 function (Review). *Mol. Med. Rep.* **20(4)**, 2979-2989 (2019).
2. Katagiri, T., Tsukamoto, S., and Kuratani, M. Accumulated knowledge of activin receptor-like kinase 2 (ALK2)/activin A receptor, type 1 (ACVR1) as a target for human disorders. *Biomedicines* **9(7)**, 736 (2021).
3. Matsuzaki, K., Xu, J., Wang, F., *et al.* A widely expressed transmembrane serine/threonine kinase that does not bind activin, inhibin, transforming growth factor  $\beta$ , or bone morphogenic factor. *J. Biol. Chem.* **268(17)**, 12719-12723 (1993).
4. Shore, E.M., Xu, M., Feldman, G.J., *et al.* A recurrent mutation in the BMP type I receptor ACVR1 causes inherited and sporadic fibrodysplasia ossificans progressiva. *Nat. Genet.* **38(5)**, 525-527 (2006).
5. Smith, K.A., Joziassse, I.C., Chocron, S., *et al.* Dominant-negative ALK2 allele associates with congenital heart defects. *Circulation* **119(24)**, 3062-3069 (2009).
6. Ensan, D., Smil, D., Zepeda-Velázquez, C.A., *et al.* Targeting ALK2: An open science approach to developing therapeutics for the treatment of diffuse intrinsic pontine glioma. *J. Med. Chem.* **63(9)**, 4978-4996 (2020).

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