

# PRODUCT INFORMATION



## TEAD1 Cofactor-Binding Domain (human, recombinant)

Item No. 35888

### Overview and Properties

**Synonyms:** NTEF-1, SV40, TEA Domain Family Member 1, TEA Domain Transcriptional Factor 1, TEF-1, Transcriptional Enhancer Factor 1

**Source:** Active recombinant human N-terminal GST- and C-terminal His-tagged TEAD1 cofactor-binding domain expressed in *E. coli*

**Amino Acids:** 209-426

**Uniprot No.:** P28347

**Molecular Weight:** 52 kDa

**Storage:** -80°C (as supplied)

**Stability:** ≥1 year

**Purity:** ≥85% estimated by SDS-PAGE

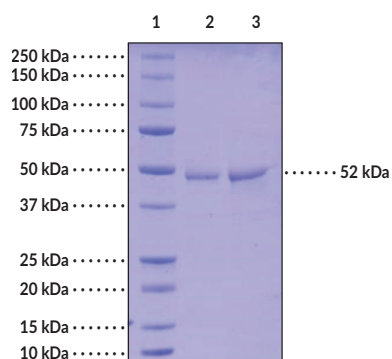
**Supplied in:** 50 mM HEPES pH 7.0, with 100 mM sodium chloride, 1 mM DTT, and 10% glycerol

**Protein**

**Concentration:** *batch specific* mg/ml

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

### Images



Lane 1: MW Markers  
Lane 2: TEAD1 Cofactor-Binding Domain (2 µg)  
Lane 3: TEAD1 Cofactor-Binding Domain (4 µg)

SDS-PAGE Analysis of TEAD1 Cofactor-Binding Domain.

Representative gel image shown; actual purity may vary between each batch.

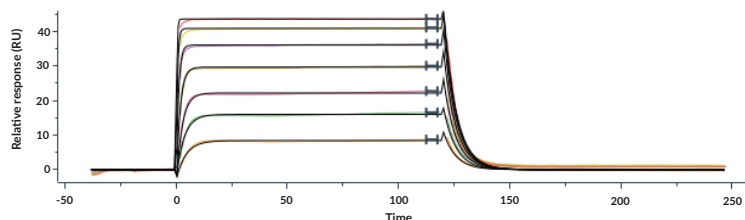


Figure 1. TEAD1 was immobilized on a Series S Sensor Chip CM5 via amine coupling. SPR analysis was used to determine YAP1 binding affinity on a Biacore™ 8K, using single-cycle kinetics with five concentrations of YAP1 (2.4-1500 nM).

**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, 09/11/2025

**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

TEA domain transcription factor 1 (TEAD1) is a transcription factor and member of the TEA domain-containing family of transcription factors.<sup>1</sup> It is composed of an N-terminal TEA DNA-binding domain, a hydrophobic proline-rich domain, and a C-terminal cofactor-binding domain (CBD), which binds TAZ, YES-associated transcriptional regulator (YAP), vestigial-like proteins (VGLLs), and several other cofactors. Lipidation of the TEAD1 CBD increases TAZ- or YAP-dependent transcriptional activity.<sup>2</sup> TEAD1 is expressed in late embryonic development and adult tissues.<sup>3</sup> It is involved in the Hippo signaling pathway, which regulates embryonic and organ development, proliferation, and cell death.<sup>1</sup> TEAD1 is specifically involved in cardiomyocyte and vascular smooth muscle cell proliferation, myogenesis, gliosis, osteoclastogenesis, and progenitor stemness, among other processes.<sup>4,5</sup> TEAD1 containing a tyrosine-to-histidine mutation at position 421 (TEAD1<sup>Y421H</sup>) is associated with Sveinsson's chorioretinal atrophy.<sup>6</sup> Increased nuclear tumor levels of TEAD1 are associated with decreased disease-free survival in patients with pancreatic ductal adenocarcinoma (PDAC).<sup>7</sup> Cayman's TEAD1 Cofactor-Binding Domain (human, recombinant) protein can be used for binding assays.

## References

1. Landin-Malt, A., Benhaddou, A., Zider, A., *et al.* An evolutionary, structural and functional overview of the mammalian TEAD1 and TEAD2 transcription factors. *Gene* **591**(1), 292-303 (2016).
2. Noritsugu, K., Suzuki, T., Dodo, K., *et al.* Lysine long-chain fatty acylation regulates the TEAD transcription factor. *Cell Rep.* **42**(4), 112388 (2023).
3. Kaneko, K.J. and DePamphilis, M.L. Regulation of gene expression at the beginning of mammalian development and the TEAD family of transcription factors. *Dev Genet.* **22**(1), 43-55 (1998).
4. Currey, L., Thor, S., and Piper, M. TEAD family transcription factors in development and disease. *Development* **148**(12), dev196675 (2021).
5. Figeac, N., Mohamed, A.D., Sun, C., *et al.* VGLL3 operates via TEAD1, TEAD3 and TEAD4 to influence myogenesis in skeletal muscle. *J. Cell Sci.* **132**(13), jcs225946 (2019).
6. Fossdal, R., Jonasson, F., Kristjansdottir, G.T., *et al.* A novel TEAD1 mutation is the causative allele in Sveinsson's chorioretinal atrophy (helicoid peripapillary chorioretinal degeneration). *Hum. Mol. Genet.* **13**(9), 975-981 (2004).
7. Drexler, R., Fahy, R., Küchler, M., *et al.* Association of subcellular localization of TEAD transcription factors with outcome and progression in pancreatic ductal adenocarcinoma. *Pancreatology* **21**(1), 170-179 (2021).

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