

# PRODUCT INFORMATION



## EED (human recombinant)

Item No. 10628

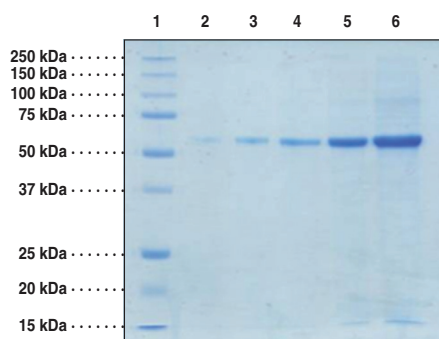
### Overview and Properties

**Synonyms:** Embryonic Ectoderm Development, WAIT-1, WD protein associating with integrin cytoplasmic tails 1  
**Source:** Recombinant N-terminal hexahistidine-tagged protein, expressed in Sf21 cells using a baculovirus expression system  
**Amino Acids:** 1-441 (full-length)  
**Uniprot No.:** O75530

Batch specific information can be found on the Certificate of Analysis or by contacting Technical Support

**Molecular Weight:** 53.5 kDa  
**Storage:** -80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein  
**Stability:** ≥9 months  
**Purity:** *batch specific* (≥95% estimated by SDS-PAGE)  
**Supplied in:** *batch specific*  
**Protein**  
**Concentration:** *batch specific*

### Image



Lane 1: MW Standards  
Lane 2: Pure EED (0.5 µg)  
Lane 3: Pure EED (1 µg)  
Lane 4: Pure EED (2 µg)  
Lane 5: Pure EED (5 µg)  
Lane 6: Pure EED (10 µg)

Representative gel image shown; actual purity may vary between each batch but protein will be ≥95% pure.

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, 05/18/2016

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

---

The human-Embryonic ectoderm development (EED) protein is a WD40 repeat-containing protein that forms part of the Polycomb repressive complex 2 (PRC2).<sup>1,2</sup> The PRC2 core complex contains the catalytic subunit EZH2 (Enhancer of zeste 2), SUZ12 (Suppressor of zeste) and EED. The EED subunit does not contain methyltransferase activity. However, transcriptional repression by PRC2-mediated trimethylation of lysine 27 on Histone H3 (H3K27me3) has been shown to be dependent on EED binding to repressive histone marks.<sup>3</sup>

## References

---

1. Schumacher, A., Lichtarge, O., Schwartz, S., *et al.* The murine polycomb-group gene *eed* and its human orthologue: Functional implications of evolutionary conservation. *Genomics* **54**, 79-88 (1998).
2. Suganuma, T. and Workman, J.L. WD40 repeats arrange histone tails for spreading of silencing. *J. Mol. Cell Biol.* **2**, 81-83 (2010).
3. Margueron, R., Justin, N., Ohno, K., *et al.* Role of the polycomb protein EED in the propagation of repressive histone marks. *Nature* **461**, 762-767 (2009).

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