
EED Protein

(Catalog No. E24030)

Description

Human Embryonic Ectoderm Development (hEED), also known as WAIT-1, GenBank Accession No. NM_003797, a.a. 2-end with N-terminal FLAG-tag.

Molecular Weight (kDa)

51 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 90 ug/ml Flag peptide, and 10% glycerol.

Source

Expressed in Sf9 cells via a baculovirus expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity>90%

Application Notes

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Ordering Information:

Products	Size	Cat. No.
EED Protein	50 µg	E24030 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.