

# PRODUCT INFORMATION



## Contactin-5 Long Isoform (human, recombinant)

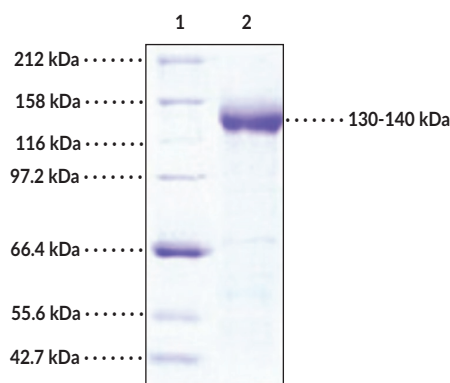
Item No. 43689

### Overview and Properties

**Synonyms:** CNTN5, hNB-2, Neural Recognition Molecule NB-2  
**Source:** Recombinant human C-terminal His-tagged contactin-5 long isoform expressed in HEK293 cells  
**Amino Acids:** 16-1,059  
**Uniprot No.:** O94779  
**Molecular Weight:** 116 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** ≥90% 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  
**Bioactivity:** See figure for details

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

### Image



Lane 1: MW Markers  
Lane 2: Contactin-5 Long Isoform

**SDS-PAGE Analysis of Contactin-5 Long Isoform.** This protein has a calculated molecular weight of 116 kDa, and a predicted N-terminus of Glu19 after signal peptide cleavage. It has an apparent molecular weight of approximately 130-140 kDa by SDS-PAGE under reducing conditions due to glycosylation.

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, 06/16/2025

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

---

Contactin-5 is a neural cell adhesion molecule and member of the immunoglobulin (Ig) superfamily.<sup>1</sup> It is composed of six Ig-like repeats, four fibronectin type III-like domains, and a hydrophobic C-terminal sequence that contains a glycosylphosphatidylinositol (GPI) anchor site. Contactin-5 is primarily expressed in the occipital lobe, amygdala, cerebral cortex, frontal lobe, thalamus, and temporal lobe of the brain but is also expressed in the thyroid.<sup>2</sup> It is involved in neurodevelopment, arborization, and sensory terminal function and interacts with amyloid- $\beta$  precursor-like protein 1 (APLP-1), protein tyrosine phosphatase  $\gamma$  (PTP $\gamma$ ), and contactin-associated protein-like 4 (CNTNAP4).<sup>1,3</sup> A SNP in *CNTN5*, the gene encoding contactin-5, is associated with increased risk and disease progression in patients with sporadic Alzheimer's disease.<sup>2</sup> Cayman's Contactin-5 Long Isoform (human, recombinant) protein consists of 1,052 amino acids, has a calculated molecular weight of 116 kDa, and a predicted N-terminus of Glu19 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 130-140 kDa due to glycosylation.

## References

---

1. Oguro-Ando, A., Zuko, A., Kleijer, K.T.E., *et al.* A current view on contactin-4, -5, and -6: Implications in neurodevelopmental disorders. *Mol. Cell. Neurosci.* **81**, 72-83 (2017).
2. Kamei, Y., Takeda, Y., Teramoto, K., *et al.* Human NB-2 of the contactin subgroup molecules: Chromosomal localization of the gene (*CNTN5*) and distinct expression pattern from other subgroup members. *Genomics* **69**(1), 113-119 (2000).
3. Shimoda, Y. and Watanabe, K. Contactins: Emerging key roles in the development and function of the nervous system. *Cell Adh. Migr.* **3**(1), 64-70 (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