

# PRODUCT INFORMATION



## CISD1 Cytosolic Domain (human, recombinant)

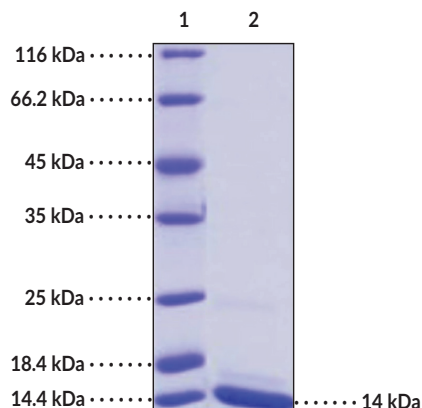
Item No. 32034

### Overview and Properties

**Synonyms:** CDGSH Iron Sulfur Domain 1, C10orf70, MitoNEET, ZCD1  
**Source:** Recombinant human N-terminal His-tagged CISD1 expressed in *E. coli*  
**Amino Acids:** 32-108  
**Molecular Weight:** 11.3 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** ≥90% estimated by SDS-PAGE  
**Supplied in:** Lyophilized from sterile 20 mM Tris, with 150 mM sodium chloride and 10% glycerol  
**Endotoxin Testing:** <1.0 EU/μg, determined by the LAL endotoxin assay

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

### Image



Lane 1: MW Markers

Lane 2: CISD1 Cytosolic Domain

#### SDS-PAGE Analysis of CISD1 Cytosolic Domain.

This protein has a calculated molecular weight of 11.3 kDa. It has an apparent molecular weight of approximately 14 kDa by SDS-PAGE under reducing conditions.

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, 06/02/2020

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

# PRODUCT INFORMATION



## Description

---

CDGSH iron sulfur domain 1 (CISD1), also known as mitoNEET, is an iron-sulfur outer mitochondrial membrane protein and member of the NEET protein family with roles in redox sensing, iron uptake, and respiratory capacity.<sup>1,2</sup> It is a homodimeric protein comprised of a  $\beta$ -cap domain and a cluster binding domain with each monomer consisting of an N-terminal in-organelle domain, a transmembrane domain, and a C-terminal cytosolic domain.<sup>1</sup> The CISD1 C-terminal domain houses the [2Fe-2S] cluster binding site which has roles in tuning redox properties of the protein in response to environmental conditions. CISD1 mRNA expression is induced by the ferroptosis inducer erastin (Item No. 17754) in HepG2 cells in an iron-dependent manner and stabilization of the CISD1 [2Fe-2S] cluster by pioglitazone (Item Nos. 71745 | 22263 | 10028) inhibits mitochondrial iron uptake, lipid peroxidation, and subsequent ferroptosis in the same cells.<sup>2</sup> Overexpression of CISD1 decreases oxidative stress-induced apoptosis in an HL-1 cardiomyocyte model of hypoxia and reoxygenation injury.<sup>3</sup> Levels of CISD1 mRNA are decreased in subcutaneous and visceral adipose tissue isolated from patients with obesity and visceral adipose tissue levels of CISD1 mRNA are positively correlated with insulin sensitivity in patients with morbid obesity.<sup>4</sup> Knockdown of *Cisd1* in mice induces mitochondrial dysfunction and increased levels of mitochondrial reactive oxygen species (ROS), as well as a loss of striatal dopamine, shortened stride length in a gait analysis, and decreased latency to fall in the rotarod test, all of which serve as markers of a Parkinsonian phenotype.<sup>5</sup> Cayman's CISD1 Cytosolic Domain (human, recombinant) protein consists of 95 amino acids and has a calculated molecular weight of 11.3 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass is approximately 14 kDa.

## References

---

1. Karmi, O., Marjault, H.-B., Pesce, L., *et al.* The unique fold and lability of the [2Fe-2S] clusters of NEET proteins mediate their key functions in health and disease. *J. Biol. Inorg. Chem.* **23(4)**, 599-612 (2018).
2. Yuan, H., Li, X., Zhang, X., *et al.* CISD1 inhibits ferroptosis by protection against mitochondrial lipid peroxidation. *Biochem. Biophys. Res. Commun.* **478(2)**, 838-844 (2016).
3. Habener, A., Chowdhury, A., Echtermeyer, F., *et al.* MitoNEET protects HL-1 cardiomyocytes from oxidative stress mediated apoptosis in an *in vitro* model of hypoxia and reoxygenation. *PLoS One* **11(5)**, e0156054 (2016).
4. Moreno-Navarrete, J.M., Moreno, M., Ortega, F., *et al.* CISD1 in association with obesity-associated dysfunctional adipogenesis in human visceral adipose tissue. *Obesity (Silver Spring)* **24(1)**, 139-147 (2016).
5. Geldenhuys, W.J., Benkovic, S.A., Lin, L., *et al.* MitoNEET (CISD1) knockout mice show signs of striatal mitochondrial dysfunction and a Parkinson's disease phenotype. *ACS Chem. Neurosci.* **8(12)**, 2759-2765 (2017).

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