

PRODUCT INFORMATION



Calreticulin (human, recombinant; His-tagged)

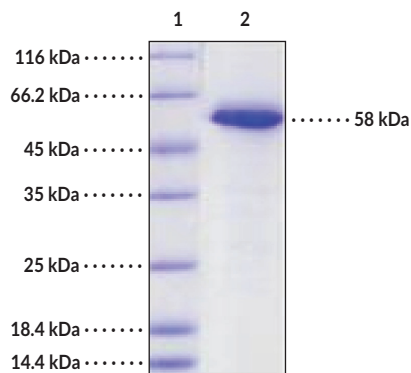
Item No. 42567

Overview and Properties

Synonyms: CAB-63, CaBP3, CALR, Calregulin, Calsequestrin-like Protein, cC1qR, CRP55, Endoplasmic Reticulum Resident Protein 60, ERp60, FLJ26680
Source: Recombinant human C-terminal His-tagged calreticulin expressed in HEK293 cells
Amino Acids: 18-413
Uniprot No.: P27797
Molecular Weight: 47.4 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥95% 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
Protein
Concentration: *batch specific* mg/ml

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

Image



Lane 1: MW Markers
Lane 2: Calreticulin

SDS-PAGE Analysis of Calreticulin. This protein has a calculated molecular weight of 47.4 kDa. It has an apparent molecular weight of approximately 58 kDa by SDS-PAGE under reducing conditions due to glycosylation and the acidity of the calreticulin C-terminal region.

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, 01/06/2025

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

Calreticulin is a calcium-binding protein involved in intracellular calcium storage and glycoprotein folding, among other processes.^{1,2} It is composed of an N-terminal globular N domain, a proline-rich P domain containing a high-affinity, low-capacity calcium-binding site, an acidic C domain containing a low-affinity, high-capacity calcium-binding site, and a C-terminal endoplasmic reticulum retention sequence.³ Calreticulin is widely expressed and localizes primarily to the lumen of the endoplasmic or sarcoplasmic reticulum in non-muscle or muscle cells, respectively, but is also found at the cell surface, in the extracellular matrix, and in the cytoplasm.^{1,3,4} It has a high calcium-binding capacity and is involved in the regulation of intracellular calcium homeostasis, which makes it an integral factor in many calcium-related processes, such as muscle contraction and the cellular stress response.^{3,4} Calreticulin also acts as a chaperone for newly synthesized glycoproteins in conjunction with protein disulfide-isomerase A3, also known as ERp57, and calnexin in the calreticulin/calnexin cycle to ensure proper glycoprotein folding.¹ Insertion or deletion mutations in exon 9 of CALR, the gene encoding calreticulin, are found in a subset of patients with essential thrombocythemia or primary myelofibrosis and are associated with increased overall survival and a lower risk of thrombosis.⁵ Cayman's Calreticulin (human, recombinant; His-tagged) protein consists of 407 amino acids, has a calculated molecular weight of 47.4 kDa, and a predicted N-terminus of Glu18 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 58 kDa due to glycosylation and the acidity of the calreticulin C-terminal region.

References

1. Wang, Q., Groenendyk, J., and Michalak, M. Glycoprotein quality control and endoplasmic reticulum stress. *Molecules* **20**(8), 13689-13704 (2015).
2. Gold, L.I., Eggleton, P., Sweetwyne, M.T., et al. Calreticulin: Non-endoplasmic reticulum functions in physiology and disease. *FASEB J.* **24**(3), 665-683 (2010).
3. Michalak, M., Milner, R.E., Burns, K., et al. Calreticulin. *Biochem J.* **285**(Pt 3), 681-692 (1992).
4. Wang, W.-A., Groenendyk, J., and Michalak, M. Calreticulin signaling in health and disease. *Int. J. Biochem. Cell Biol.* **44**(6), 842-846 (2012).
5. Klampfl, T., Gisslinger, H., Harutyunyan, A.S., et al. Somatic mutations of calreticulin in myeloproliferative neoplasms. *N. Engl. J. Med.* **369**(25), 2379-2390 (2013).

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