

PRODUCT INFORMATION



Mcl-1L Cytoplasmic Domain (human, recombinant)

Item No. 33983

Overview and Properties

Synonyms: Bcl-2-like Protein 3, EAT/Mcl-1, Induced Myeloid Leukemia Cell Differentiation Protein
Source: Active recombinant human N-terminal His-tagged MCL1 expressed in *E. coli*
Amino Acids: 2-327
Uniprot No.: Q07820
Molecular Weight: 36 kDa
Storage: -80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
Stability: ≥6 months
Purity: ≥90% estimated by SDS-PAGE
Supplied in: 40 mM Tris-HCl, pH 8.0, with 110 mM sodium chloride, 2.2 mM potassium chloride, 0.04% Tween 20, 3 mM DTT, 20% glycerol, and 200 mM imidazole

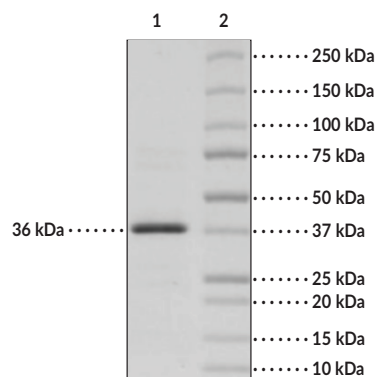
Protein

Concentration: *batch specific* mg/ml

Activity: See figure for details

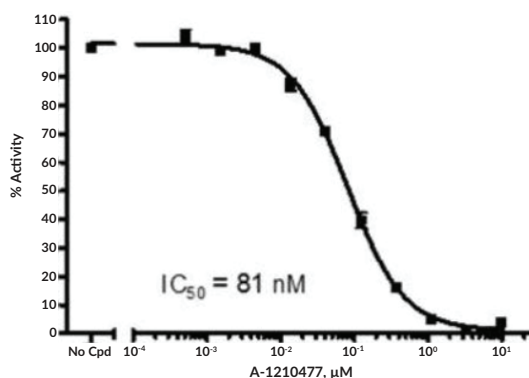
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: Mcl-1L Cytoplasmic Domain
Lane 2: MW Markers

SDS-PAGE Analysis of Mcl-1L Cytoplasmic Domain.



Inhibition of MCL-1 by A-1210477 (Item No. 21113).

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, 11/22/2021

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

Myeloid cell leukemia-1 (Mcl-1) is a Bcl-2 family member protein that is involved in the regulation of apoptosis.¹ Alternative splicing of *MCL1* pre-mRNA produces three isoforms, anti-apoptotic Mcl-1 long (Mcl-1L) and pro-apoptotic Mcl-1 short and Mcl-1 extra short (Mcl-1S and Mcl-1ES, respectively). Mcl-1L is composed of a C-terminal helix that acts as a membrane anchor, four Bcl-2 homology (BH) domains, and a PEST domain. The BH1, BH2, and BH3 domains facilitate heterodimerization with pro-apoptotic proteins, including Bax and Bak, and homodimerization with Mcl-1S and Mcl-1ES.¹⁻³ Mcl-1L is expressed during embryogenesis and hematopoiesis, as well as during nervous and immune system development, and localizes to the mitochondrial membrane.^{4,5} It inhibits apoptosis by preventing Bax and Bak oligomerization, which form a pore on the outer mitochondrial membrane that allows release of cytochrome c into the cytoplasm.¹ Knockdown of *MCL1* sensitizes oral cancer cells to cisplatin and tumor levels of Mcl-1L are increased relative to adjacent normal tissues in patients with oral cancer.⁶ Overexpression of *MCL1* is associated with poor prognosis in several cancers, including solid tumors and hematological malignancies.¹ Cayman's Mcl-1L Cytoplasmic Domain (human, recombinant) protein can be used for enzyme activity assays.

References

1. Wang, H., Guo, M., Wei, H., *et al.* Targeting MCL-1 in cancer: Current status and perspectives. *J. Hematol. Oncol.* **14**(1), 67 (2021).
2. Bae, J., Leo, C.P., Hsu, S.Y., *et al.* MCL-1S, a splicing variant of the antiapoptotic BCL-2 family member MCL-1, encodes a proapoptotic protein possessing only the BH3 domain. *J. Biol. Chem.* **275**(33), 25255-25261 (2000).
3. Kim, J.-H., Sim, S.-H., Ha, H.-J., *et al.* MCL-1ES, a novel variant of MCL-1, associates with MCL-1L and induces mitochondrial cell death. *FEBS Lett.* **583**(17), 2758-2764 (2009).
4. Opferman, J.T. and Kothari, A. Anti-apoptotic BCL-2 family members in development. *Cell Death Differ.* **25**(1), 37-45 (2018).
5. Yang, T., Kozopas, K.M., and Craig, R.W. The intracellular distribution and pattern of expression of Mcl-1 overlap with, but are not identical to, those of Bcl-2. *J. Cell Biol.* **128**(6), 1173-1184 (1995).
6. Palve, V., Mallick, S., Ghaisas, G., *et al.* Overexpression of Mcl-1L splice variant is associated with poor prognosis and chemoresistance in oral cancers. *PLoS One* **9**(11), e111927 (2014).

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