

Cofilin 1 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # C37-54H

Lot # D061-4

Product Description

Recombinant full-length human Cofilin 1 was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_005507](#).

Gene Aliases

CFL1, CFL

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 2mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

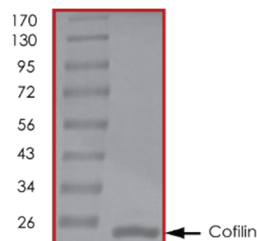
Scientific Background

Cofilin 1 (non-muscle), also known as CFL1, is a widely distributed intracellular actin-modulating protein that binds and depolymerizes filamentous F-actin and inhibits the polymerization of monomeric G-actin in a pH-dependent manner (1). It is involved in the translocation of actin-cofilin complex from cytoplasm to nucleus (2). CFL1 processed proteins are most probably produced by HIV-1 protease digestion.

References

1. Maciver, S.K. et al: The ADF/cofilin family: actin-remodeling proteins." *Genome Biol.* 2002; 3 (5): reviews 3007.
2. Samstag, Y. Et al: Interaction of cofilin with the serine phosphatases PP1 and PP2A in normal and neoplastic human T lymphocytes." *Adv. Enzyme Regul.* 2004;97-211.

Purity



The purity of Cofilin 1 was determined to be **>95%** by densitometry.
Approx. MW **23kDa**.

Cofilin 1 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog #	C37-54H
Lot #	D061-4
Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.