

Recombinant Human CD40 Ligand (CD40LG)

Recombinant human protein expressed in E. coli cells

Catalog # C810-40N

Lot # G3211-16

Product Description

Recombinant Human CD40 Ligand (CD40LG) was expressed in E. coli cells. The protein accession number is [P29965](#)

Gene Aliases

Soluble CD40 Ligand, Tumor necrosis factor ligand superfamily member 5, TNFSF5, TNF-related activation protein, TRAP, CD154, gp39, T-BAM, CD40, T cell antigen Gp39, CD-40L

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

Lyophilized from a 0.2 µm filtered solution in PBS (pH 7.0)

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

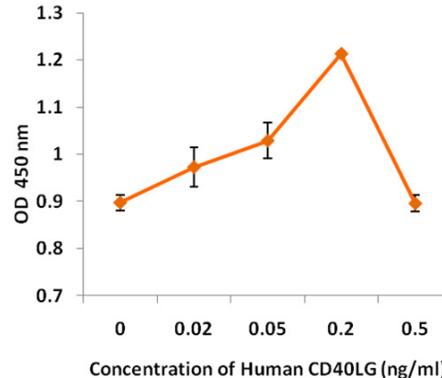
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

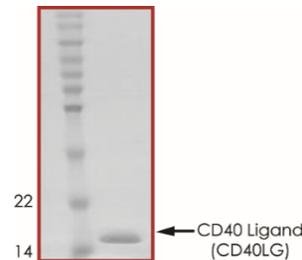
CD40LG protein is expressed on all B lymphocytes during various stages of development, activated T cells and monocytes, follicular dendritic cells, thymic epithelial cells, and various carcinoma cell lines. It is expressed on most mature B cell malignancies and on some early B cell acute lymphocytic leukemias. Human CD40LG has been found to be the receptor for a protein involved in an X-linked immunodeficiency syndrome. Based on its homology with other members of the TNF ligand protein superfamily, CD40LG is also referred to as TNFSF5. The soluble form of CD40LG is generated by the intracellular proteolytic processing of the full length CD40LG. Recombinant Human CD40LG is a 17 kDa protein that represents the soluble segment of full length CD40LG and encompasses the receptor binding TNF-like domain.

Specific Activity



The ED50 as determined by dose-dependent proliferation of acute myeloid leukemia was found to be <5 ng/ml

Purity



Recombinant Human CD40 Ligand (CD40LG) resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **17.0 kDa**

Recombinant Human CD40 Ligand (CD40LG)

Recombinant human protein expressed in E. coli cells

Catalog Number	C810-40N
Specific Lot Number	G3211-16
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

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.