

Catalog #	Aliquot Size
U866-40N-15	15 µg
U866-40N-100	100 µg
U866-40N-1000	1 mg

Recombinant Human Uteroglobin (SCGB1A1)

Recombinant human protein expressed in E. coli cells

Catalog # U866-40N

Lot # G3223-24

Product Description

Recombinant Human Uteroglobin (SCGB1A1) was expressed in E. coli cells. The protein accession number is [P11684](#)

Gene Aliases

Clara cell phospholipid-binding protein, CCPBP, Clara cells 10 kDa secretory protein, CC10, Secretoglobin family 1A member 1, Urinary protein 1, Urine protein 1, UP1

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

Reconstitution Protocol

Reconstitute at 100 µg/mL in sterile PBS

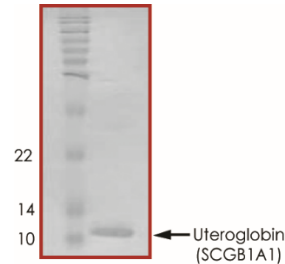
Storage and Stability

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

UP1 is a non-glycosylated secretoglobin family member protein which induces anti-inflammatory and anti-tumorigenic responses. The expression of Uteroglobin is induced by steroid hormones such as estrogen which is further enhanced by the non-steroid hormone prolactin. UP1 has been shown to induce growth stimulatory effects, sequester pro-inflammatory mediators and carcinogens and inhibit cell migration, cell invasion, platelet aggregation and T cell differentiation.

Purity



Recombinant Human Uteroglobin (SCGB1A1) resolved on a 17% SDS_PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **8.2 kDa**

Recombinant Human Uteroglobin (SCGB1A1)

Recombinant human protein expressed in E. coli cells

Catalog Number	U866-40N
Specific Lot Number	G3223-24
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.