

| Catalog # | Aliquot Size |
|----------------------|---------------|
| T813-40N-20 | 20 µg |
| T813-40N-100 | 100 µg |
| T813-40N-1000 | 1 mg |

Recombinant Human TFF3

Recombinant human protein expressed in E. coli cells

Catalog # T813-40N

Lot # G3223-15

Product Description

Recombinant Human TFF3 was expressed in E. coli cells. The protein accession number is [Q07654](#)

Gene Aliases

Intestinal trefoil factor, ITF, Polypeptide P1.B, hP1.B
Recombinant Human Trefoil Factor 3 (TFF3)

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.

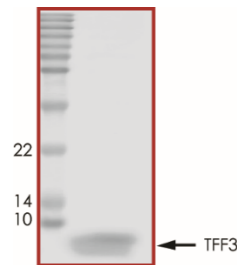
Scientific Background

Found in the gastrointestinal tract, the three Trefoil proteins, TFF1, 2, and 3 have been shown to play an important role in intestinal mucosal defence and repair. In addition to being expressed by goblet cells and in the uterus, TFF3 has also been shown to express in certain cancers providing a potential use as a molecular marker for those cancers.

Specific Activity

Determined by its ability to chemoattract human MCF-7 cells using a concentration range of 1.0-10.0 µg/ml

Purity



Recombinant Human TFF3 resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **6.8 kDa**

Recombinant Human TFF3

Recombinant human protein expressed in E. coli cells

| | |
|---------------------|---|
| Catalog Number | T813-40N |
| Specific Lot Number | G3223-15 |
| 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.