

TARC (human), (recombinant)

Thymus and activation regulated chemokine (TARC), also known as CCL17, is a chemokine produced by thymus tissue constitutively and activated PBMCs (mainly DCs). TARC signals through the CCR4 receptor to induce chemotaxis of Th2 cells. TARC is thought to be important in asthma and allergic diseases, along with bacterial and viral infections.

Manuals, SDS & CofA

[View Online »](#)

- High purity
- Carrier-free

Handling & Storage

Use/Stability	It is recommended that a carrier protein (0.1% HSA or BSA) is added for long term storage.
Handling	Centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	CCL17, Thymus and activation regulated chemokine, ABCD-2, SCYA17
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the ability to chemoattract human T cells which is typically observed at a concentration of 2-40ng/ml.
Endotoxin Content	≤1 EU/μg protein measured by kinetic LAL analysis.
Formulation	Lyophilized from 0.02% TFA. Sterile filtered.
MW	~8.1kDa
Purity	≥98% (HPLC, Reducing and Non-reducing SDS-PAGE, UV spectroscopy at 280 nm)
Reconstitution	When reconstituting the product, gently pipet and wash down the sides of the vial to ensure full recovery of the protein into solution. It is recommended to reconstitute the lyophilized product with sterile water at a concentration of 0.1 mg/ml, which can be further diluted into other aqueous solutions.
Source	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 71 amino acids.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com