

Chemerin monoclonal antibody (Tigy-1)

Chemerin (tazarotene-induced gene 2; TIG2) is a chemotactic agent that was identified as the ligand for the orphan G protein-coupled receptor ChemR23. Chemerin has been shown to be expressed abundantly in adipose tissue and to function as an adipokine. Chemerin potentiates insulin-stimulated glucose uptake and regulates adipogenesis and adipocyte metabolism.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-868-C100	100µg
------------------	-------

Manuals, SDS & CoFA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	RARRES2, Retinoic acid receptor responder protein 2, TIG2, Tazarotene-induced gene 2 protein, RAR-responsive protein
Application	ELISA, WB
Clone	Tigy-1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant rat chemerin (aa 23-157).
Isotype	IgG2a
Purity Detail	Protein G-affinity purified.
Recommendation	Western blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Dilutions/Conditions	
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human, Rat



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

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#).

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.