

# 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 a 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

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

**Regulatory Status** RUO - Research Use Only

## Product Details

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

UniProt ID

Q5BK77

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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