

# Calcitonin gene related peptide polyclonal antibody

Calcitonin gene-related peptide (CGRP) is a 37-amino acid disulphide-bridged sensory neuropeptide present in the central and peripheral nervous systems of many animal species. Two forms of CGRP ( $\alpha$  and  $\beta$ ) have been identified from both rat and human, and the sequence of chicken CGRP has also been reported, showing striking structural similarity to the mammalian forms. CGRP is abundant in perivascular sensory neurons and is known to be a potent vasodilator. Recent work supports the rôle of CGRP in mediating skin vasodilatation and differential mesenteric artery vasorelaxation, as well as serving as an endogenous myocardial protective substance with a possible involvement in the inhibition of human platelet aggregation. CGRP is frequently co-localised with Substance P and immunoelectron microscopic observations have shown that the peptides are co-stored in single secretory granules both in sensory neurons and in thyroid C-cells. There are many hundreds of reports in the scientific literature describing the localisation and co-existence of CGRP with other transmitter molecules in a vast array of mammalian and non-mammalian species.

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

Citations: 29

[View Online »](#)

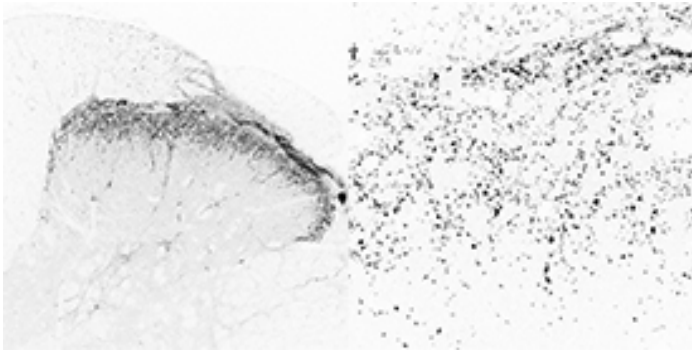
## Ordering Information

[Order Online »](#)

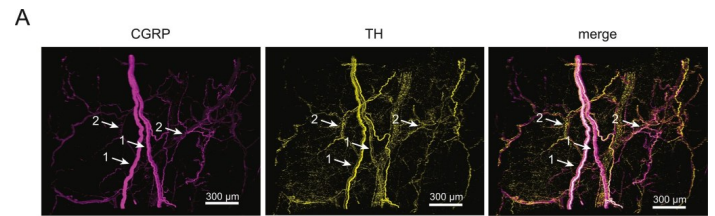
BML-CA1137-0025	25µl
BML-CA1137-0100	100µl

Manuals, SDS & CofA

[View Online »](#)

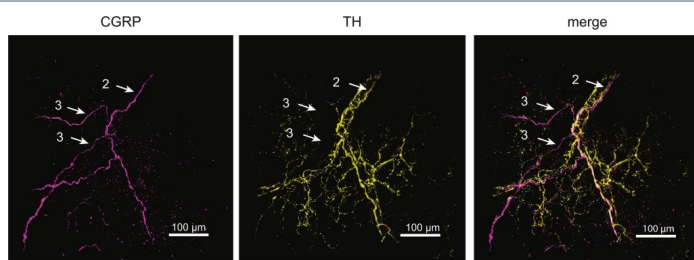


Immunohistochemistry analysis: Vibratome (70µm) sections of rat spinal cord immunostained with sheep polyclonal antiserum to CGRP (Prod. No. BML-CA1137) at a dilution of 1:1000. Negative images of fluorescence detection procedure (Alexa488 conjugated secondary antibody). Micrograph courtesy of Professor Andrew Todd (University of Glasgow).



(A) 2D representative of a 3D reconstruction of an inguinal WAT (iWAT) section immunostained with calcitonin gene-related peptide (CGRP, magenta) and the tyrosine hydroxylase (TH, yellow) that illustrates how a large nerve bundle subdivides into smaller nerve bundles which in turn penetrate the tissue alongside blood vessels. Nerve bundle (1), innervation along blood vessel (2). Scale bar= 300 µm. (B) 2D representative of a 3D reconstruction of the iWAT parenchyma of both CGRP-positive (magenta) and the TH-positive (yellow) neuronal networks in iWAT. Innervation along blood vessel (2), parenchymal innervation (3). Scale bar= 100 µm.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Adipose mTORC2 is essential for sensory innervation in white adipose tissue and whole-body energy homeostasis. *Mol Metab* (2022)



(A) 2D representative of a 3D reconstruction of an inguinal WAT (iWAT) section immunostained with calcitonin gene-related peptide (CGRP, magenta) and the tyrosine hydroxylase (TH, yellow) that illustrates how a large nerve bundle subdivides into smaller nerve bundles which in turn penetrate the tissue alongside blood vessels. Nerve bundle (1), innervation along blood vessel (2). Scale bar= 300 µm. (B) 2D representative of a 3D reconstruction of the iWAT parenchyma of both CGRP-positive (magenta) and the TH-positive (yellow) neuronal networks in iWAT. Innervation along blood vessel (2), parenchymal innervation (3). Scale bar= 100 µm.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Adipose mTORC2 is essential for sensory innervation in white adipose tissue and whole-body energy homeostasis. *Mol Metab* (2022)

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	CGRP
Application	IHC
Crossreactivity	Likely to react with other mammalian species (based on homology, not tested).
Formulation	Liquid. In PBS containing 0.1% sodium azide.
Host	Sheep
Immunogen	Synthetic peptide corresponding to a portion of rat $\alpha$ -calcitonin gene-related peptide (CGRP).
Purity Detail	Ammonium sulphate precipitation.
Source	Purified from sheep serum.
Species Reactivity	Human, Mouse, Rabbit, Rat
Technical Info / Product Notes	<p>Application Data details.</p> <p>Test tissues: Rat thoracolumbar spinal cord.</p> <p>Fixatives: Recommended fixative is 4% formaldehyde (prepared from para-polymer) in 0.1M sodium phosphate buffer (pH 7.2-7.4). Perfusion fixation is recommended if possible.</p>
UniProt ID	P01256

### Worry-free Guarantee

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



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)