Calcitonin gene related peptide monoclonal antibody (CD8)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-CA1135-0100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CGRP

Application IHC

Clone CD8

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of C-terminal

rat calcitonin gene-related peptide (CGRP).

lgG1

Purity Detail Partially purified.

Species Reactivity Guinea pig, Human

UniProt ID P01256

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

