

CGRP Affinity SorbENT

ORDERING INFORMATION

Catalogue Number: A19482

Stability: ≥ 2 years

Storage: +4°C

Shipping: ice pack

**For research laboratory
use only.**

**Not for human
diagnostic use.**

PRODUCT DESCRIPTION

The following is the purification protocol for CGRP sorbent provided by SPI-BIO.

- A. Pour the gel in an empty column (around 10 x 1 cm).
- B. Let the storage solution elute entirely (the gel should become dry but not for a long time).
- C. Add around 20 mL of plasma on top of the column with the outlet of the column closed.
- D. Let it incubate overnight at +4°C with gentle agitation.
- E. Elute the plasma that is now free of CGRP and ready to be used to dilute standards and samples in the assay.
- F. So as to regenerate the gel (eliminate the bound CGRP), wash the column with around 20 mL of 0.1 M HCl pH 1.5 and then with around 50 mL of 0.1 M Phosphate buffer pH 7.4.
- G. You can then store the sorbent in 3 to 4 mL of EIA buffer (buffer of the kit) at +4°C till the next use. It is stable at +4°C for many months.

Related Products

1. Human CGRP EIA kit ref. #A05481.
2. Rat CGRP EIA kit ref. #A05482.

Should you wish further information, do not hesitate to contact us.



• Parc d'Activités du Pas du Lac • 10 bis avenue Ampère - F-78180 Montigny Le Bretonneux Cedex
Tel: 33 (0)1 39 30 62 60 • Fax: 33 (0)1 39 30 62 99 • E-Mail: sales@spibio.com
Web : www.spibio.com

