

Unblocked protein A beads

Cat. No.	Format
C03020004	880 µl
C03020003	2.8 ml
C03020002	8 ml

Description

The unblocked protein A beads are agarose beads that have been extensively validated in chromatin immunoprecipitation assays (ChIP). These beads are intended for isolation of immune complexes (chromatin and specific antibody) in ChIP experiments. The beads are suitable for immunoprecipitation of rabbit polyclonal Abs, mouse IgG2a, IgG2b and IgA, guinea pig IgG, dog IgG, pig IgG. The beads should be washed before use.

Format

Supplied as a suspension in PBS with 20% of ethanol. Keep the beads in liquid suspension during storage as drying will result in reduced performance.

Storage

Store at 4°C. Do not freeze. Stable for 1 year.

Precautions

This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: July, 2023