

DNA precipitant

ChIP and DNA IP grade reagent

Separately available

Cat. No. C03030002

Size: 1000 µl

Concentration: 3 M solution (Supplied as a 3 M stock solution (pH 5.2). 50 µl allow to precipitate IP'd DNA from a volume of 500 µl)

Storage buffer: PBS containing 0.5 mM PMSF and 50% glycerol.

Storage: The product is stable for at least 2 years when stored at 4°C.

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

Description

This IP grade product is used for the precipitation of DNA after chromatin immunoprecipitation (ChIP) and after methylated DNA immunoprecipitation (Methyl DNA IP). It is included in our **Transcription ChIP KIT**, in our **Histone ChIP KIT** and in our **Methylated DNA Immunoprecipitation KIT - MeDIP**.

It is now separately available. Following IP and washes, the immunoprecipitated (IP'd) material is submitted to Phenol/Chloroform/Isomanyl alcohol (PCI) extraction. Then, to the aqueous phase obtained after PCI extraction of the IP'd material, our DNA precipitant is added with **IP grade DNA co-precipitant** and ice-cold 96% ethanol. For more details about these kits, see instruction manual on our web site (www.diagenode.com).

Diagenode sa. BELGIUM | EUROPE

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

Diagenode Inc. 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: 10.2007