

## 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](http://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](mailto:orders@diagenode.com)  
[info@diagenode.com](mailto: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](mailto:orders.na@diagenode.com)  
[info.na@diagenode.com](mailto:info.na@diagenode.com)

Last update: 10.2007