

PRODUCT NAME		
PCR-grade water		
ChIP grade reagent Separately available		
Cat. No: kch-508-028	Size: 28 ml	Lot #: D005
Cat. No: kch-508-085	Size: 85 µl	Lot #: D005

Description: PCR-grade water is designed for use in all molecular biology applications. It is 0.1 µm membrane-filtered for DNase and RNase activity.

This product is used to dilute the purified DNA from the Immunoprecipitated (IP'd) chromatin. It is included in our **OneDay ChIP KIT (cat# kch-onedIP-060, -180)** and our **LowCell# ChIP KIT (cat# kch-maglow-016)**. It is now separately available. For more details about these kits, see instruction manual on our web site (www.diagenode.com).

Format: Supplied as a solution ready to use.

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.

Availability date: October 29, 2007

Last data sheet update: October 26, 2007

Lot #: D005/ aliquoting day: October 02, 2007