

PRODUCT NAME		
PCR-grade water		
ChIP grade reagent Separately available		
Cat. No: <b>kch-508-028</b>	Size: 28 ml	Lot #: D005
Cat. No: <b>kch-508-085</b>	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](http://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