

PRODUCT NAME			
Proteinase K			
ChIP grade reagent Separately available			
Cat. No: kch-507-100	Size: 100 µl	Concentration: 100 x stock solution	Lot #: D004
Cat. No: kch-507-250	Size: 250 µl	Concentration: 100 x stock solution	Lot #: D004

Description: Proteinase K is a non-specific serine protease. This ChIP grade product has been extensively validated to isolate the 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 100 x stock solution.

Storage: The product is stable for at least 2 years when stored at -20°C. Avoid multiple freeze-thaw cycles.

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 #: D004/ aliquoting day: October 02, 2007