

| PRODUCT NAME                               |              |                                     |             |
|--|--------------|-------------------------------------|-------------|
| Proteinase K                               |              |                                     |             |
| ChIP grade reagent<br>Separately available |              |                                     |             |
| Cat. No: <b>kch-507-100</b>                | Size: 100 µl | Concentration: 100 x stock solution | Lot #: D004 |
| Cat. No: <b>kch-507-250</b>                | 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](http://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