

<b>PRODUCT NAME</b> DNA co-precipitant		
ChIP and DNA IP grade reagent Separately available		
Cat. No: <b>kch-216-100</b>	Size: 100 µl	Concentration: 2 µg/µl stock solution

**Description:** This IP grade product is optimized to be used as carrier for DNA purification after chromatin immunoprecipitation (ChIP) and after methylated DNA immunoprecipitation (Methyl DNA IP). It is included in our **Transcription ChIP KIT (cat# kch-redTBP-012)**, in our **Histone ChIP KIT (cat# kch-orgHIS-012)** and in our **Methylated DNA Immunoprecipitation KIT-MeDIP kit (cat# mc-green-003)**.

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 DNA material, our DNA co-precipitant is added with **IP grade DNA precipitant (cat# kch-217-001)** 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)).

**Format:** Supplied as a 2 µg/µl stock solution. 10 µg allow to precipitate IP'd DNA from a volume of 500 µl.

**Storage:** The product is stable for 1-2 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 #:** D003/ aliquoting day: October 02, 2007