

PRODUCT NAME			
GenDNA RNase (DNase-free)			
Methyl DNA IP grade reagent Separately available			
Cat. No: mcg-605-010	Size: 10 µl	Concentration: 500 x stock solution	Lot #: D001

Description: Our RNase is DNase-free. No boiling treatment is needed; it can be used directly from the vial. This product is particularly well suited for the use in methylated DNA isolation procedures. It is included in our **XL GenDNA Extraction MODULE (cat# mc-magme-003)** and it is now separately available. For more details about this kit, see instruction manual on our web site (www.diagenode.com).

Format: 500 x stock solution ready to use. Residual RNA can be removed by adding 2 µl of GenDNA RNase (DNase-free) per ml of DNA sample and incubating for 1 hour at 37°C.

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 #: D001/ aliquoting day: October 02, 2007