

## CATS paired-end sequencing primer

**Cat. No.** C17011050-50/500

**Format:** 50 µl/500 µl

**Product description:**

This primer is specifically designed to be used in paired-end sequencing with CATS Library Preparation on Illumina platforms (MiSeq, HiSeq).

**Format:** This product contains a custom primer for paired-end sequencing at a concentration of 100 µM in 50 µl or 500 µl of ultrapure water.

**Storage:** Store at -20°C. Avoid multiple freeze/thaw cycle.

**Shipment:** Dry ice.

**Shelf-life:** 2 years.

**Precautions:** This product is for research use only. Not for use in diagnostics or therapeutic procedures.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** August 26, 2016