

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 RNA-seq libraries on Illumina platforms (MiSeq, HiSeq).

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