

## Mouse SFI1 promoter primer pair

Cat. No. **C17021046**

Product Name	Quantity	Concentration	Storage
Cat. No. C17021046-50	50 µl	5 µM each	-20°C/-4°F

### Product description

The primer pair Cat. No. C17021046 is specific to SFI1 Centrin Binding Protein from mouse. SFI1 is a protein coding gene that has been identified as being hydroxymethylated using DNA immunoprecipitation assays (hMeDIP). The primers are optimized to be used in quantitative polymerase chain reaction (qPCR).

### Applications

*For capture of hydroxymethylated DNA (e.g. hMeDIP - Hydroxymethylated DNA Immunoprecipitation)*

The region amplified with SFI1 primer pair corresponds to a hydroxymethylated region in mouse. The primers can be used as a positive control to monitor the success of the capture by qPCR. The size of the expected amplicon is 86 bp.

**Storage conditions:** Store at -20°C/-4°F. For long storage, store at -80°C /-112°F. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic 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: December 16, 2020