

## Bisulfite conversion reagent for RRBS

Cat. No. **C02030035**

Format: 8 rxns

### Content

Description	Quantity	Storage
BS Conversion Reagent	1 tube	Room temperature
BS Dilution Buffer	300 µl	Room temperature
BS Solubilization Buffer	790 µl	Room temperature
BS Reaction Buffer	160 µl	Room temperature

### Product description

This reagent is compatible with the Premium RRBS Kit and provides reagents for additional bisulfite conversion reactions. Please check the Premium RRBS Kit manual for further information.

Reduced Representation Bisulfite Sequencing provides a powerful method to efficiently analyze DNA methylation at the **single nucleotide level** without the higher costs associated with whole genome bisulfite sequencing. By cutting the genome using the restriction MspI enzyme (CCGG target sites) followed by size selection, DNA is enriched to represent CpG-rich regions (including CpG islands), in which DNA methylation marks are typically found. Thus, RRBS provides a **cost-effective** method for analyzing DNA methylation by reducing the part of the genome that actually needs to be sequenced and focusing on relevant CpG islands.

#### 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: 01.2017