

## MethylTaq Plus 2X Master Mix

**Cat. No.** C09010012

**Format:** 135 µL

**Shipping conditions:** Ship on dry ice or ice packs, depending on the country of destination.

**Storage conditions:** Store at -15°C to -25°C.

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

### Description

MethylTaq Plus 2X Master Mix is a ready-to-use cocktail containing all components required for the amplification of bisulfite-converted NGS libraries, except primers and template. This convenient 2X Master Mix contains MethylTaq Plus DNA polymerase, each dNTP (dATP, dCTP, dGTP, dTTP, but not dUTP) [0.3 mM], MgCl<sub>2</sub> [2.5 mM] and reaction buffer.

MethylTaq Plus DNA polymerase is a high-performance Hot Start thermostable recombinant DNA polymerase designed for the PCR amplification of uracil-containing templates with high efficiency, low bias and high fidelity. This highly robust enzyme produces excellent results in demanding applications and it is recommended for PCR after bisulfite conversion. MethylTaq Plus DNA polymerase requires a 5 minutes initial denaturation step at 95°C to reach maximal initial activity. Because of the composition of the reaction buffer, a denaturation temperature of 98°C is recommended during the cycles. Optimal cycle numbers should be determined empirically.

#### 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:** September 1, 2020