

Data Sheet

MECP2

Human recombinant, with N-terminal GST-tag
Catalog #: 50250
Lot#: 91028 **Conc.:** 2.35 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
 124 mM NaCl, 2.2 mM KCl, 18 mM
 glutathione, 10% glycerol and 3 mM DTT.

Stability: >6 months at -80 °C

References:

1. Amir, R.E., *et al.* (1999). *Nat. Genet.* **23** (2): 185–8.
2. Lewis, J.D., *et al.* (1992). *Cell* **69** (6): 905–14.
3. Chahrour M, *et al.* (2008). *Science* **320** (5880): 1224–9.

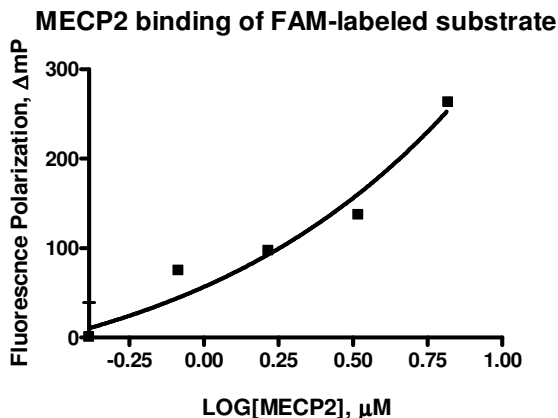
Description:

Human recombinant Methyl CpG binding Protein 2 (Rett syndrome) (MECP2), GenBank Accession # NM_004992), a.a. 78-162 with N-terminal GST-tag, MW= 36 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of DNA binding, screening inhibitors, and selectivity profiling.

Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 10 μg MECP2
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 36 kDa
Purity: ≥90%

