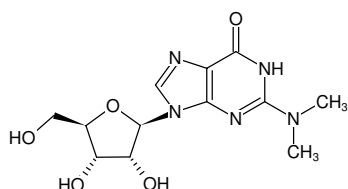




## N<sup>2</sup>,N<sup>2</sup>-Dimethylguanosine

(m<sup>2</sup>,2G)

Cat. No.	Amount
N-1087-10	10 mg
N-1087-100	100 mg



Structural formula of N<sup>2</sup>,N<sup>2</sup>-Dimethylguanosine

### For research use only!

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 4 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 48 months after date of delivery

**Molecular Formula:** C<sub>12</sub>H<sub>17</sub>N<sub>5</sub>O<sub>5</sub>

**Molecular Weight:** 311.29 g/mol

**CAS#:** 2140-67-2

**Purity:** ≥ 98 % (HPLC)

**Form:** white to off-white solid

**Spectroscopic Properties:** λ<sub>max</sub> 262 nm; ε 12.3 L mol<sup>-1</sup> cm<sup>-1</sup> (pH 11.0)

### Selected References:

Dai *et al.* (2017) Selective Enzymatic Demethylation of N<sup>2</sup>,N<sup>2</sup>-Dimethylguanosine in RNA and Its Application in High-Throughput tRNA Sequencing. *Angew. Chem. Int. Ed. Engl.* **56** (18):5017.

Su *et al.* (2014) Quantitative analysis of ribonucleoside modifications in tRNA by HPLC-coupled mass spectrometry. *Nat. Protoc.* **9** (4):828.

Edqvist *et al.* (1995) Enzymatic formation of N<sup>2</sup>,N<sup>2</sup>-dimethylguanosine in eukaryotic tRNA: importance of the tRNA architecture. *Biochimie* **77** (1):54.