

## Poly T gripNA™ Probe

---

**Catalog No:** 29007, 29008

### **Quality Control:**

Poly T gripNA™ Probe has been optimized to ensure efficient recovery of mRNA from cells, tissue and Total RNA. It is a hetero-oligomer constructed of alternating thymine (T) pPNA (phosphono PNA) and thymine (T) HypNA (hydroxy PNA) monomers in a 1:1 ratio with a biotin molecule at the 5' end. pPNA monomers contain N-(2-hydroxyethyl)phosphonoglycine or N-(2-aminoethyl)phosphonoglycine as the backbone, while HypNA monomers contain trans-4-hydroxy-L-proline as the backbone molecule.

Poly T gripNA Probe is included in Active Motif's mTRAP™ mRNA Isolation Kits and is suitable for use in various poly A<sup>+</sup> mRNA isolation procedures.

Poly T gripNA Probe is guaranteed to be at least 70% biotinylated. 4-5 nmol of Poly T gripNA Probe is sufficient to isolate up to 150 µg of mRNA from either cell or tissue lysates. 600 pmol is sufficient to isolate up to 10 µg of mRNA from 500 µg of Total RNA.

### **Storage:**

Poly T gripNA Probe is supplied lyophilized. It should be stored at 4°C for short-term and -20°C for long-term storage. Catalog No. 29007 should be resuspended in 500 µl of the supplied HPLC water. Catalog No. 29008 should be resuspended in 400 µl of the supplied HPLC water. After resuspension, the storage conditions remain the same as those for the lyophilized probe.