

## pAM Forward primer

---

**Catalog No:** 53049

**Format:** 250 pmol

**Quality Control:** Primers are purified by gel electrophoresis or HPLC.

### Background

pAM Forward and Reverse primers are designed for use with the pAM vector system by Active Motif. These primers can be used for DNA sequencing to confirm cloning reactions with the pAM\_1C empty vector (Catalog No. 53023).

pAM Forward: 5'-GTTCCGAAAGTTGCCTTTTA-3'

pAM Reverse: 5'-CGGATCCTTATCGATTTTACCAC-3'

### Contents

pAM Forward and Reverse primers are provided lyophilized. Reconstitute with sterile distilled water or TE Buffer (pH 7.5-8.0). Once reconstituted, store at -20°C.

**Storage and Guarantee:** Lyophilized product shipped at RT. Store at -20°C

This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.