

TE, pH 8.0

Catalog No: 37515

Format: 35 ml

Quality Control: Active Motif's TE, pH 8.0 (part number 104890) is a commonly used buffer solution in molecular biology. TE buffer can be used for a variety of applications including solubility of DNA or RNA solutions. TE buffer is also used to wash Protein G agarose beads prior to use in immunoprecipitation.

TE buffer is 10 mM Tris-HCl, 1 mM EDTA, pH 8.0 and is provided ready to use. If using TE to prepare Protein G agarose beads (Catalog No. 37499), we recommend using an equal volume of TE, pH 8.0 to the volume of Protein G agarose beads. Invert to mix and centrifuge at 1250 x g for 1 minute. Remove the supernatant volume equivalent to the amount of TE added. Repeat wash steps if desired. Beads are ready to use in ChIP, Co-IP or other applications.

Storage and Guarantee: Store at room temperature. 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.