

mTRAP™ Lysis Buffer

Catalog No: 29011

Quality Control:

Composition

200 mM Tris, pH 7.5	200 mM NaCl
500 mM GTC	Proprietary mix of detergents

Active Motif's Lysis Buffer is supplied in the mTRAP Kits. It consists of the chaotropic reagent guanidine thiocyanate (GTC) and a mixture of detergents that insure the complete inactivation of RNases in cell and tissue lysates. The composition of mTRAP Lysis Buffer has been optimized for use with Active Motif's Poly T PNA Probe (Catalog Nos. 29007 & 29008) for isolation of poly A⁺ mRNA.

Active Motif's Lysis Buffer is suitable for lysing any type of cell or tissue, while maintaining RNA integrity. It is suitable for tissues high in RNases (e.g. placenta tissue) and tissues that may require vigorous mechanical homogenization, such as plant tissues.

Additional Materials Required:

Protease (Catalog No. 29011) or some similar proteinase (e.g. Proteinase K) should be added to the Lysis Buffer to a final concentration of 200-400 µg/ml prior to use. The final concentration of Protease may need to be optimized for each type of cell lysate.

Use 15 ml Lysis Buffer for up to 1 g of tissue or 2 x 10⁸ mammalian cells. For complete information, obtain mTRAP product manuals from one of our Technical Service departments, or download one at www.activemotif.com.

Storage:

Store at room temperature. If a white precipitate develops, heat the buffer at 37°C for 10 minutes prior to use.

This product is guaranteed for 6 months from date of receipt.

Caution

Lysis Buffer contains the corrosive guanidine thiocyanate, a strong denaturant. Exercise appropriate caution when working with this solution, including the use of gloves, a labcoat and protective eyewear.