

Lysis Buffer AM1

Catalog No: 37512

Format: 50 ml

Quality Control: Active Motif's Lysis Buffer AM1 (part number 100566) is used to prepare nuclear and whole-cell lysates from mammalian cell culture or tissue samples. Lysis Buffer AM1 contains phosphatase (but not protease) inhibitors.

Lysis Buffer AM1 is provided as a 1X solution. We recommend supplementing the buffer with Protease Inhibitor Cocktail (Catalog No. 37490) prior to use. Lysis Buffer AM1 is compatible with Active Motif's Nuclear Extract Kit (Catalog No. 40010) and TransAM Transcription Factor-DNA Binding assays.

Storage and Guarantee: Lysis Buffer AM1 contains phosphatase inhibitors which gives it a yellow coloration. The product is shipped at room temperature and should be stored at 4°C upon arrival. 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.