

Extraction Buffer

Catalog No: 37513

Format: 100 ml

Quality Control: Active Motif's Extraction Buffer (part number 102168) is used for the isolation of core histone proteins from cell culture and tissue samples. The highly basic nature of histone proteins often precludes them from isolation using traditional methods for protein extraction. Instead, histones must be isolated under acidic conditions to retain solubility.

Extraction Buffer is provided as a ready to use solution. We recommend incubating a homogenous cell solution in ice-cold Extraction Buffer (1.5 ml per 150 mm dish) for 2 hours to overnight at 4°C. Following extraction, centrifuge the cells at maximum speed for 5 minutes and transfer the supernatant to a new tube. Add 1/4 volume of 5X Neutralization Buffer (Catalog No. 37514) to the supernatant (*e.g.* add 0.25 ml 5X Neutralization Buffer to 1 ml crude histones for a final volume of 1.25 ml neutralized histones). This provides a crude core histone extraction.

For purified histone extraction, we suggest using Active Motif's Histone Purification Kits (Catalog Nos. 40025, 40026 & 40027).

Storage and Guarantee: Extraction Buffer is shipped at room temperature and can 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.