

5X Neutralization Buffer

Catalog No: 37514

Format: 50 ml

Quality Control: Active Motif's 5X Neutralization Buffer (part number 102169) is used in 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.

5X Neutralization Buffer is provided as a 5X stock solution. We recommend using it in combination with Extraction Buffer (Catalog No. 37513) to prepare crude histone lysates. Incubate 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: 5X Neutralization 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.