

## Core Histones (Chicken)

**CATALOG NO.:** HMT-35-435

**LOT NO.:**

**DESCRIPTION:** Core histones (including histone H5) purified from chicken erythrocytes by a modification of the method of Schnitzler<sup>1</sup> followed by acid extraction/dialysis<sup>2</sup>.

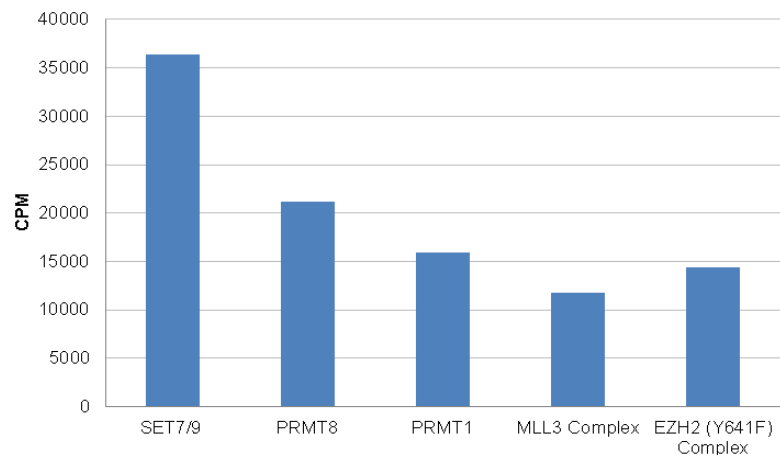
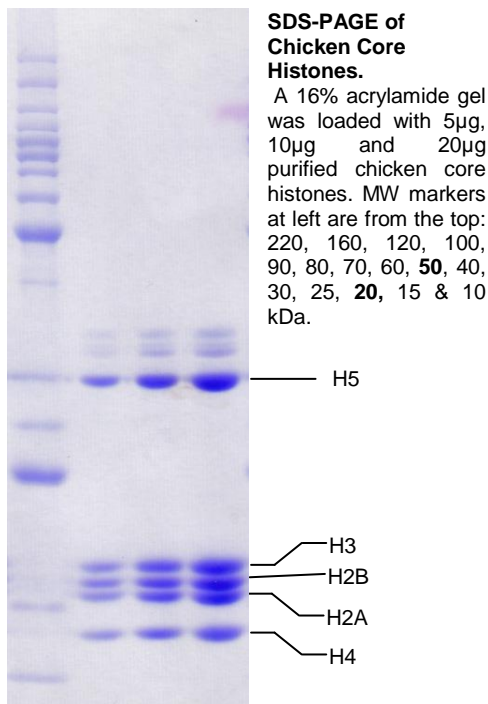
**PURITY:** >90% by SDS-PAGE.

**APPLICATIONS:** Useful for the assay of various histone methyltransferases (see figure below) by methods employing radiolabeling with [<sup>3</sup>H]-S-adenosylmethionine (SAM) (e.g. gel electrophoresis/autoradiography or filterplate/scintillation counting). Standard HMT reaction conditions: 50 mM Tris-HCl, pH 8.5, 50 mM NaCl, 5 mM MgCl<sub>2</sub> with 0.05 mg/mL core histones and 1 μM [<sup>3</sup>H]-SAM as substrates.

**SUPPLIED AS:** \_\_\_ μg/μl in dH<sub>2</sub>O as determined by Bradford protein assay.

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted portion should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 μl) or storage of diluted solutions is not recommended.

**REFERENCE:** 1) G. Schnitzler *Current Protocols in Molecular Biology* 2000 21.5.1-21.5.12, 2) D. Shechter, *et al Nature Protocols* 2007 (2)6: 1445-1457.



**Histone Methyltransferase Activity on Chicken Core Histones.** Assays (25 μL) were performed with a scintillation/filter plate assay. Incubations were 60 min., 30°C with 20nM SET7/9 (RBC Cat. # HMT-11-133); 100nM PRMT8 (RBC Cat. # HMT-11-135), 75nM PRMT1 (RBC Cat. #HMT-11-119), 200nM MLL3 Complex (RBC Cat. #HMT-15-107), or 50nM EZH2 (Y641F) Complex (RBC Cat. #HMT-25-173) plus 0.05mg/mL chicken core histones and 1 μM [<sup>3</sup>H]-SAM.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.