hBSA: Cationized with Hexamethylenediamine

Catalog Number:5601Size:10 mgMolecular Weight:67 kD

Solubility: Soluble in water

Description: hBSA is prepared by treating the native BSA with hexamethylenediamines that replace most of negatively-

charged carboxyl groups with positively-charged primary amines, resulting a highly positively-charged hBSA. The cationization significantly increases the immunogenicity compared to native BSA. In addition, the increased number of primary amines provides more conjugation sites available for hapten molecules with general conjugation methods. The modification of BSA with hexamethylenediamine provides a longer space

between the carrier protein and the hapten.

Storage/Stability: Store at -20°C/1 year

Format: Lyophilized in PBS, pH 7.2.

Immunogen: Use as a carrier protein for immunization

Purification: BSA is purified by a fractionation method, and is supplied with the purity over 97% by SDS.

Applications: hBSA itself acts an excellent immnogen with a greater immunogenicity compared to the native BSA. With

increased number of free amines, more antigen molecules can be coupled to hBSA. When a stronger immunogenicity and a high concentration of hapten are needed, hBSA is a good choice for the immunogen preparation of small hapten molecules, particularly for a longer space between the carrier protein and the

hapten.

References: Sheng-Liang Deng, Ping Li, Hong-Bin Liu, Shu-Ming Yang (2014) Preparation and characterization of

ultrasensitive and specific polyclonal antiserum against ciprofloxacin based on cationized bovine serum

albumin. Chemical Papers 68 (11) 1505-1513.

Technical Support: support@aatbio.com; Tel: 408-733-1055