

Avanti's Water Soluble KLA with Huzzah[®]

Avanti's new HSA lipid delivery system, Huzzah[®], improves the cellular delivery of lipid species through their conjugation with human serum albumin (HSA), a physiologically relevant carrier protein. This conjugate system eliminates the need for organic solvents, which are well-known to adversely affect viability and the phenotypic characteristics of cells. Avanti's latest conjugate, Huzzah[®] KLA, provides the researcher with the ability to deliver this new synthetic glycan in aqueous solution.

KLA has been shown to be 10x more effective than lipopolysaccharide



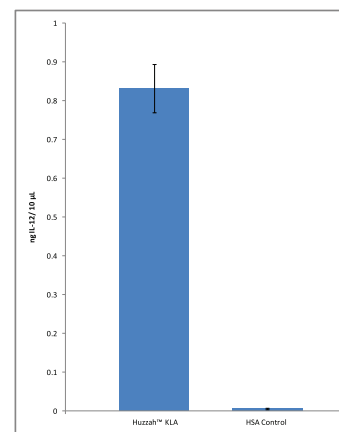
Huzzah[®] is shipped as a lyophilized powder. Photograph on right shows complex dissolved in water.

Now in Stock:

- Huzzah[®] KLA Avanti Number 360500
- Huzzah[®] S1P Avanti Number 360492
- Huzzah[®] LPA Avanti Number 360130
- Huzzah[®] NBD-So Avanti Number 360205

Adjuvant preparation without sonication or organic solvents

Many Lipid A-based adjuvant systems require lengthy sonication and serial dilution. Huzzah[®] KLA is provided as a freeze-dried conjugate with human serum albumin that only requires the addition of saline to be ready for use. Interleukin 12 (IL-12) secretion from serum-deprived J774A mouse monocyte/macrophage is shown in adjacent Figure after 24h exposure to Huzzah[®] KLA. IL-12 content of 10 μ L samples was determined by sandwich ELISA with an anti-mouse IL-12/IL-23 p40 capture antibody kit (R&D Systems, MP400) against standards of mouse IL-12 (11.7-750 pg/mL). The IL-12 concentration was determined by measuring absorbances at 450 nm and 570 nm and subtracting the A570 from the A450 values.



More details and recent data available at avantilipids.com