

BSA ELISA Kit

Cat.No. L00976

Key Features



Ultra Sensitivity

LoB: 0.14 ng/mL; LoQ: 0.5 ng/mL



High Specificity

No HSA interference

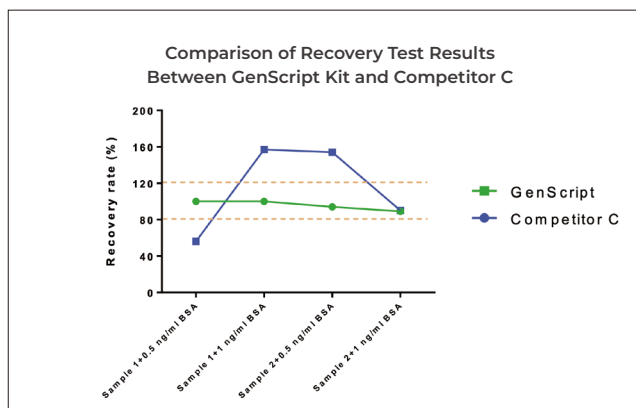


Reliable Standard

Traceable standard & accurate calibration

Competitive Performance

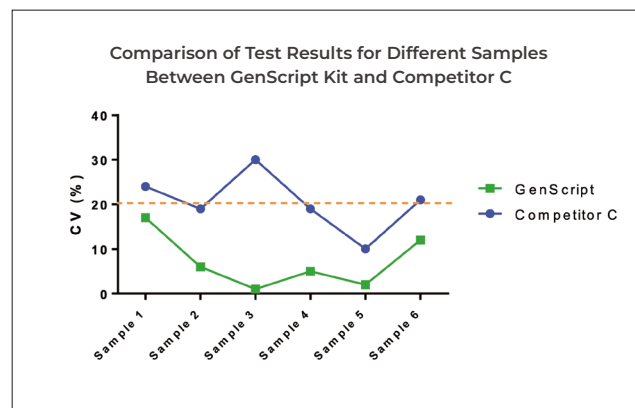
Consistently High Recovery



*Two samples were spiked with low concentration of BSA and then tested using the GenScript kit and competitor C kit to evaluate the recovery rate.

BSA ELISA kit demonstrates a recovery rate of 80%–120%, ensuring greater accuracy.

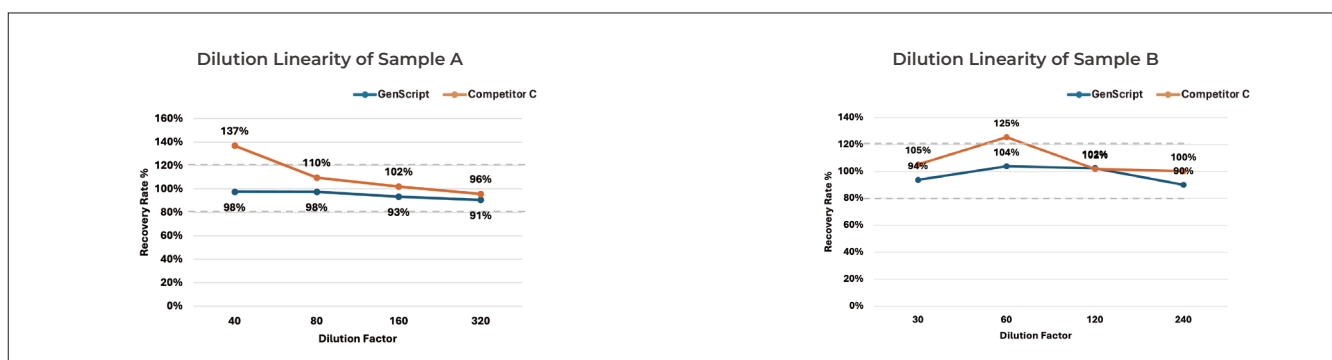
Low Variation



*Different samples were tested using GenScript kit and competitor C kit to evaluate the coefficient of variation (CV%).

The test results from the GenScript kit exhibit a CV of less than 10%, ensuring more stable outcomes.

Wide Dilution Linearity

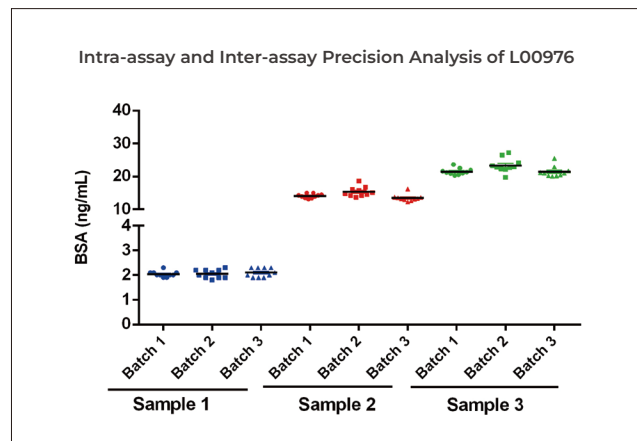


*GenScript BSA ELISA kit and competitor C kit were used to test two identical samples at different dilutions to evaluate the recovery rate.

The results from the GenScript kit remained consistent and unaffected by the dilution factor.

Validation Data

• High Precision



Intra-assay and Inter-assay CV $\leq$ 10%.

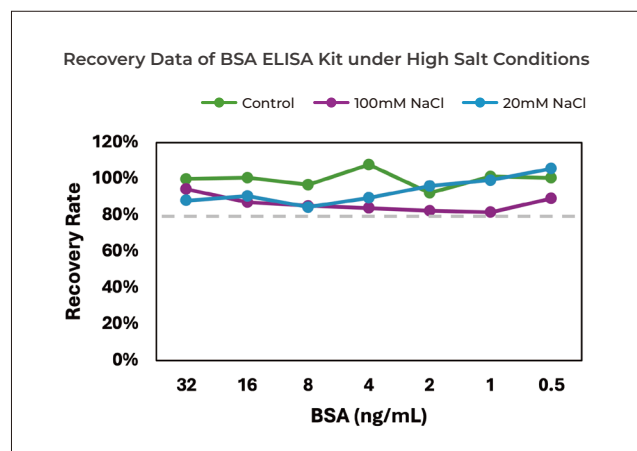
• High Specificity

HSA Spiked Conc. (ng/mL)	BSA Theoretical Conc. (ng/mL)	BSA Measured Conc. (ng/mL)	BSA Recovery Rate (%)
20	8	6.5	81
10	8	6.8	83
1	8	7.7	96
0.1	8	8.5	107

*Different concentrations of HSA (Human Serum Albumin) were spiked into the samples and blank matrices to evaluate the specificity of L00976.

The results of L00976 are not affected by the HSA.

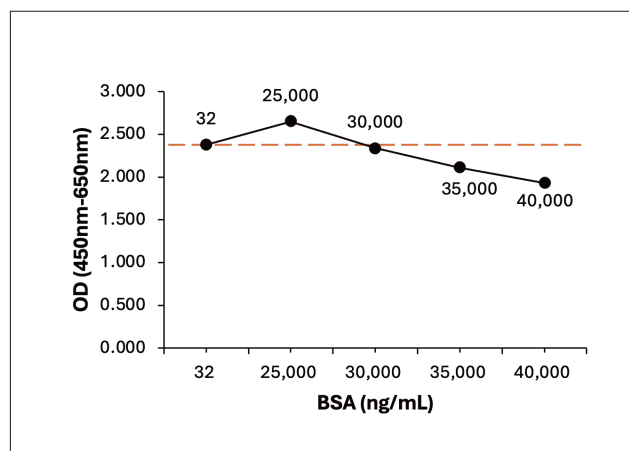
• High Salt Tolerance



*BSA ELISA Kit was used to test samples diluted with 100mM NaCl or 20mM NaCl.

The results of BSA ELISA Kit are not affected by high salt condition.

• Hook Effect



The HOOK effect appears when the BSA concentration in the samples over 30,000 ng/mL.