

BSA ELISA Detection Kit

Overview

Cat.No. HG-BS001

This kit is designed for the quantitative detection of residual BSA content in intermediates, semi-finished products and finished products of various biological products by using a double-antibody sandwich method.

Specification

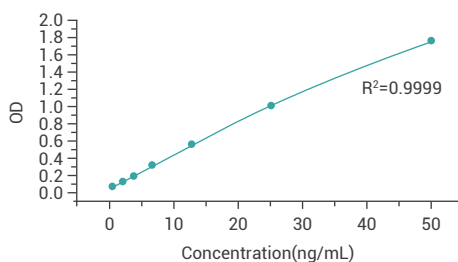
Assay range: 1.56-50 ng/mL

Limit of quantitation: 1.56 ng/mL

Limit of detection: 0.5ng/mL

Precision: CV%≤10%, RE%≤±15%

Standard curve



Datasheet

Concentration of standard (ng/mL)	OD value1	OD value2	Mean value
50	1.775	1.803	1.789
25	1.091	0.995	1.043
12.5	0.571	0.591	0.581
6.25	0.324	0.308	0.316
3.125	0.193	0.179	0.186
1.56	0.132	0.127	0.130
0	0.072	0.088	0.080