

DNase I ELISA Detection Kit

Overview

Cat.No. HG-DI001

This kit is designed for the quantitative detection of residual Dnase I content added in RNA pharmaceutical processes by using a double-antibody sandwich method.

Specification

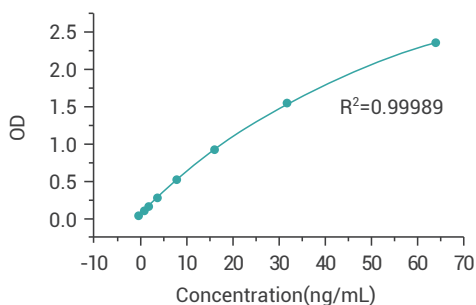
Assay range: 1-64 ng/mL

Limit of quantitation: 1 ng/mL

Limit of detection: 0.5 ng/mL

Precision: CV%≤10%, RE%≤±15%

Standard curve



Datasheet

Concentration of standard (ng/mL)	OD value1	OD value2	Mean value
64	2.373	2.347	2.360
32	1.520	1.546	1.533
16	0.933	0.934	0.9335
8	0.518	0.536	0.527
4	0.259	0.291	0.275
2	0.166	0.168	0.167
1	0.115	0.109	0.112
0	0.057	0.049	0.053