

RNase Inhibitor ELISA Detection Kit

Overview

Cat.No. HG-RI001

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

Specification

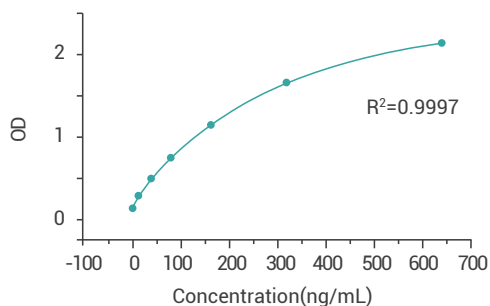
Assay range: 20-640 ng/mL

Limit of quantitation: 20 ng/mL

Limit of detection: 5 ng/mL

Precision: CV% \leq 10%, RE% \leq \pm 15%

Standard curve



Datasheet

Concentration of standard (ng/mL)	OD value1	OD value2	Mean value
640	2.213	2.110	2.162
320	1.661	1.641	1.651
160	1.181	1.143	1.162
80	0.743	0.755	0.749
40	0.486	0.478	0.482
20	0.298	0.292	0.295
0	0.151	0.15	0.151