

# Benzonase Nuclease ELISA Detection Kit

## Overview

Cat.No. HG-BE001

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

## Specification

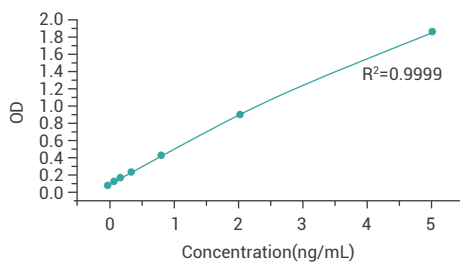
**Assay range:** 0.05-5ng/mL

**Limit of quantitation:** 0.05ng/mL

**Limit of detection:** 0.05ng/mL

**Precision:** CV% $\leq$ 10%, RE% $\leq$ ±15%

## Standard curve



## Datasheet

Concentration of standard (ng/mL)	OD value1	OD value2	Mean value
5	1.925	1.774	1.849
2	0.917	0.865	0.891
0.8	0.447	0.441	0.444
0.32	0.228	0.209	0.218
0.128	0.146	0.137	0.141
0.0512	0.114	0.107	0.110
0	0.090	0.091	0.091