

# T7 RNA Polymerase ELISA Detection Kit

## Overview

Cat.No. HG-TP001

This kit is designed for the quantitative detection of residual T7 RNA Polymerase content added in RNA pharmaceuticals processes by using a double-antibody sandwich method.

## Specification

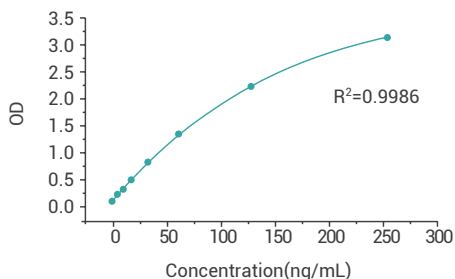
**Assay range:** 4-256 ng/mL

**Limit of quantitation:** 4 ng/mL

**Limit of detection:** 2 ng/mL

**Precision:** CV%≤10%, RE%≤±15%

## Standard curve



## Datasheet

Concentration of standard (ng/mL)	OD value1	OD value2	Mean value
256	3.109	3.183	3.146
128	2.281	2.213	2.247
64	1.411	1.394	1.4025
32	0.787	0.845	0.816
16	0.463	0.447	0.455
8	0.299	0.307	0.303
4	0.235	0.235	0.235
0	0.141	0.133	0.137