

# dsRNA ELISA Detection Kit

distributed in the US/Canada by:  
EAGLE BIOSCIENCES, INC.  
20A NW Blvd, Suite 112 Nashua, NH 03063  
Phone: 617-419-2019 FAX: 617-419-1110  
[www.EagleBio.com](http://www.EagleBio.com) info@eaglebio.com



## Overview

Cat.No. HG-DS001

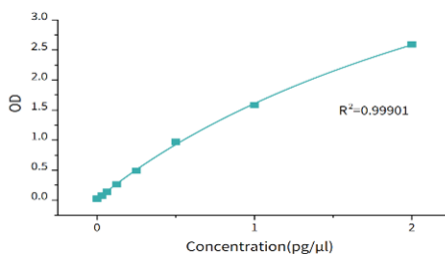
This dsRNA ELISA detection kits of BlueKit® series use a double antibody sandwich method to quantitatively detect the content of double stranded RNA (dsRNA) in samples. The detected dsRNA is 60 bp or more in length, and is not related to its nucleic acid sequence.

## Specification

Standard type	Assay range (pg/μL)	Limit of quantitation (pg/μL)	Limit of detection (pg/μL)	Recovery rate	CV%
No modification	0.0156-0.5	0.0156	0.001	80-120%	≤ 10%
N1-Me-pUTP modified	0.0312-1	0.0312	0.001	80-120%	≤ 10%
pUTP modified	0.0156-0.5	0.0156	0.001	80-120%	≤ 10%
5-OMe-UTP modified	0.0625-1	0.0625	0.01	80-120%	≤ 10%

\*Range of linearity: Coefficient of determination > 0.99

## Standard curve



## Datasheet

Concentration of standard (pg/μL)	OD value1	OD value2	Mean value
2	2.8412	2.7362	2.7887
1	1.8725	1.9135	1.8930
0.5	1.0863	1.1207	1.1035
0.25	0.623	0.6055	0.6143
0.125	0.3388	0.3292	0.3340
0.0625	0.1947	0.1885	0.1916
0.0312	0.1192	0.1247	0.1220
0	0.0567	0.0518	0.0543