

# Product Information & ELISA Manual

Equine Triiodothyronine ELISA Kit  
(Colorimetric)  
NBP2-60175

Enzyme-linked Immunosorbent Assay  
for quantitative detection.

## Contact

Web: [www.bio-techne.com/brands/novus-biologicals/](http://www.bio-techne.com/brands/novus-biologicals/)  
Email: [nb-customerservice@bio-techne.com](mailto:nb-customerservice@bio-techne.com)  
P: 888.506.6887 // P: 303.730.1950 // F: 303.730.1966

Novus kits are  
guaranteed for 6 months  
from date of receipt.

**For research use only.  
Not for diagnostic or  
therapeutic procedures.**

## PRECISION

### **Intra-assay Precision (Precision within an assay): CV%<15%**

Three samples of known concentration were tested twenty times on one plate to assess.

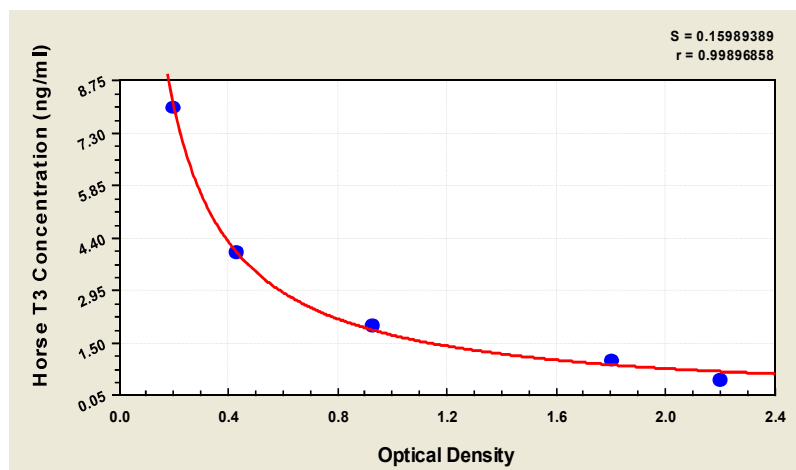
### **Inter-assay Precision (Precision between assays): CV%<15%**

Three samples of known concentration were tested in twenty assays to assess.

Sample	Intra-Assay Precision			Inter-Assay Precision		
	1	2	3	1	2	3
n	20	20	20	20	20	20
Mean(ng/ml)	1.955	1.964	1.945	1.976	1.950	2.003
SD	0.025	0.019	0.028	0.036	0.036	0.033
CV(%)	2.733	2.119	3.052	3.997	4.011	3.685

## TYPICAL DATA

These standard curves are provided for demonstration only. A standard curve should be generated for each set of samples assayed.



ng/ml	OD1	OD2	Average
0.5	2.198	2.183	2.191
1.0	1.799	1.807	1.803
2.0	0.936	0.958	0.947
4.0	0.467	0.448	0.458
8.0	0.239	0.227	0.233

## LOD

0.25ng/ml