

Product Information & ELISA Manual

Human B7-H3/CD276 ELISA Kit (Colorimetric)
NBP3-43452

Enzyme-linked Immunosorbent Assay
for quantitative detection.

Contact

Web: www.bio-techne.com/brands/novus-biologicals/
Email: 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.**

Table of Contents

1. Intended Use	3
2. Introduction	3
3. General References	4
4. Assay Principle	5
5. Handling & Storage	5
6. Kit Components	5
7. Materials Required but <i>Not</i> Supplied	6
8. General ELISA Protocol	7
8.1. Preparation and Storage of Reagents	7
8.2. Sample collection, storage and dilution	8
8.3. Assay Procedure (Checklist)	9
9. Calculation of Results	10
10. Typical Data	10
11. Performance Characteristics	11-12
12. Technical Hints and Limitations	13
13. Troubleshooting	14
14. Notes	15

1. Intended Use

The Human B7-H3/CD276 ELISA Kit (Colorimetric) is to be used for the *in vitro* quantitative determination of human B7-H3/CD276 in cell culture supernatants, serum and plasma. This ELISA Kit is for research use only.

2. Introduction

B7 family of immunoregulatory proteins is composed of ten members: B7-1 (CD80), B7-2 (CD86), B7-H1 (PD-L1), B7-DC (PD-L2), B7-H2, B7-H3, B7-H4, B7-H5 (VISTA), B7-H6 and B7-H7. B7-H3 (or CD276) is a 316aa long type I transmembrane protein. B7-H3 shares 20-27% amino acid identity with other B7 family ligands. Glycosylated B7-H3 protein has a molecular weight of approximately 100 kDa (1). A B7-H3 'soluble' isoform has been detected in plasma and B7-H3 is also expressed on exosomes (2).

B7-H3 is ubiquitously expressed by cells in the non-hematopoietic compartment, such as fibroblasts and epithelial cells, it can be induced on T cells and NK cells. Although B7-H3 expression is elevated in tumors, B7-H3 is also constitutively expressed at higher levels in the liver than in other healthy tissue (3). B7-H3 mRNA is expressed in most normal tissues, but the B7-H3 protein expression is limited in normal tissues, because of its posttranscriptional regulation by miRNAs (4). Specifically, miRNA-29 suppresses B7-H3 expression in normal tissues by targeting the B7-H3 3'-untranslated mRNA region.

Little is known about the receptors for these B7-family ligands and their downstream signaling. TLT2 (triggering receptor expressed on myeloid cells-like transcript 2) has been characterized as a putative receptor for B7-H3, but it is dispensable for T cell response, suggesting that other unknown receptors should bind B7-H3 (5). Although receptors remain unidentified, soluble B7-H3 has been shown to bind to CD4⁺ T, CD8⁺ T, NK and NKT cells and the extent of its binding is increased upon T cell activation.

B7-H3 inhibits the activation and function of T cells, potently suppressing the proliferation, cytokine production and cytotoxicity of activated T cells. It also inhibits natural killer (NK) cell activation and has a proinflammatory role leading to cytokine release from monocytes and/or macrophages. However, B7-H3 was initially characterized as a co-stimulatory molecule required for optimal promotion of T cell proliferation and cytokine production, which makes the exact role of B7-H3 controversial.

The **B7-H3/CD276 (human) ELISA Kit** detects the soluble human B7- H3 (CD276) protein that is ectopically expressed in various cancers (6) and its levels in serum of patients with cancer suggests it can be used as a non-invasive biomarker for diagnosis, prognosis and/or treatment response.

3. General References

1. B7-H3 in Cancer - Beyond Immune Regulation: F. Flem-Karlsen, et al.; Trends Cancer **6**, 401 (2018)
2. B7-H3 in Medulloblastoma-Derived Exosomes; A Novel Tumorigenic Role: I.J. Purvis, et al.; Int. J. Mol. Sci. **19**, 7050 (2020)
3. Potential Therapeutic Targets of B7 Family in Colorectal Cancer: C. Wang C. et al.; Front. Immunol. **11**, 681 (2020)
4. Regulation of cancer immune escape: The roles of miRNAs in immune checkpoint proteins: Q. Yang, et al.; Cancer Lett. **431**, 73 (2018)
5. Triggering receptor expressed on myeloid cell-like transcript 2 (TLT-2) is a counter-receptor for B7-H3 and enhances T cell responses: M. Hashiguchi, et al.; PNAS **105**, 10495 (2008)
6. The Role of CD276 in Cancers: S. Liu, et al.; Front. Oncol. **11**, 654684 (2021)

4. Assay Principle

This assay is a sandwich Enzyme Linked-Immunosorbent Assay (ELISA) for quantitative determination of human B7-H3/CD276 in cell culture supernatants, serum and plasma. A monoclonal antibody specific for human B7-H3/CD276 has been precoated onto the 96-well microtiter plate. Standards (STD) and samples are pipetted into the wells for binding to the coated antibody. After extensive washing to remove unbound compounds, B7-H3/CD276 is recognized by the addition of a monoclonal biotinylated antibody specific for human B7-H3/CD276 (DET). After removal of excess biotinylated monoclonal antibody, streptavidin-peroxidase (STREP-HRP) is added. Following a final washing, peroxidase activity is quantified using the substrate 3,3',5,5'-tetramethylbenzidine (TMB). The intensity of the color reaction is measured at 450nm after acidification and is directly proportional to the concentration of B7-H3/CD276 in the samples.

5. Handling & Storage

- Reagent must be stored at 2-8°C when not in use
- Plate and reagents should be at room temperature before use.
- Do not expose reagents to temperatures greater than 25°C.

6. Kit Components

- | | | |
|---|----------------------|--------------------|
| • 1 vial human B7-H3 Standard (lyophilized) | (100 ng) | (STD) |
| • 1 vial B7-H3/CD276 Detection Antibody | (20 µl) | (DET) |
| • 1 vial HRP Labeled Streptavidin (lyophilized) | (2 µg) | (STREP-HRP) |
| • 2 bottles Wash Buffer 10X | (2 x 30 ml) | (Wash Buffer 10X) |
| • 1 bottle ELISA Buffer 10X | (1 x 30 ml) | (ELISA Buffer 10X) |
| • 1 bottle TMB Substrate Solution | (12 ml) | (TMB) |
| • 1 bottle Stop Solution | (12 ml) | (STOP) |
| • 1 plate coated with B7-H3/CD276 Antibody | (6 x 16-well strips) | |
| • 2 plate Covers (plastic film) | | |
| • 2 silica Gel Minibags | | |

7. Materials Required but *Not* Supplied

- Microtiterplate reader at 450nm
- Calibrated precision pipettes. Disposable pipette tips
- Deionized water
- Microtubes or equivalent for preparing dilutions
- Disposable plastic containers for preparing working buffers
- Plate washer: automated or manual
- Glass or plastic tubes for diluting and aliquoting standard

8. General ELISA Protocol

8.1. Preparation and Storage of Reagents

NOTE: Prepare just the appropriate amount of the buffers necessary for the assay.

- **Wash Buffer 10X** has to be diluted with deionized water 1:10 before use (e.g. 30 ml Wash Buffer 10X + 270 ml water) to obtain Wash Buffer 1X.
- **ELISA Buffer 10X** has to be diluted with deionized water 1:10 before use (e.g. 10 ml ELISA Buffer 10X + 90 ml water) to obtain ELISA Buffer 1X.
- **Detection Antibody (DET)** has to be diluted to 1:1'000 in ELISA Buffer 1X (10 µl DET + 10 ml ELISA Buffer 1X).

NOTE: The diluted Detection Antibody is not stable and cannot be stored!

- **HRP Labeled Streptavidin (STREP-HRP)** has to be reconstituted with 100 µl of ELISA Buffer 1X.
 - After reconstitution of STREP-HRP, prepare aliquots and store them at -20°C. **Avoid freeze/thaw cycles.**
 - Dilute the reconstituted STREP-HRP to the working concentration by adding 50 µl in 10 ml of ELISA Buffer 1X (1:200).

NOTE: The diluted STREP-HRP is not stable and cannot be stored!

- **Human B7-H3/CD276 (STD)** has to be reconstituted with 100 µl of ELISA Buffer 1X.
 - This reconstitution produces a stock solution of 1 µg/ml. Mix the standard to ensure complete reconstitution and allow the standard to sit for a minimum of 15 minutes **at room temperature**. Mix well prior to making dilutions.

NOTE: The reconstituted standard is aliquoted and stored at -20°C!

- Dilute the standard protein concentrate (STD) (**1 µg/ml**) in ELISA Buffer 1X. A seven-point standard curve using 2-fold serial dilutions in ELISA Buffer 1X is recommended.
- Suggested standard points are:
20, 10, 5, 2.5, 125, 0.625, 0.3125 and 0 ng/ml.

Dilute further for the standard curve:

To obtain	Add	Into
20 ng/ml	20 µl of B7-H3/CD276 STD (1 µg/ml)	980 µl of ELISA Buffer 1X
10 ng/ml	300 µl of B7-H3/CD276 (20 ng/ml)	300 µl of ELISA Buffer 1X
5 ng/ml	300 µl of B7-H3/CD276 (10 ng/ml)	300 µl of ELISA Buffer 1X
2.5 ng/ml	300 µl of B7-H3/CD276 (5 ng/ml)	300 µl of ELISA Buffer 1X
1.25 ng/ml	300 µl of B7-H3/CD276 (2.5 ng/ml)	300 µl of ELISA Buffer 1X
0.625 ng/ml	300 µl of B7-H3/CD276 (1.25 ng/ml)	300 µl of ELISA Buffer 1X
0.03125 ng/ml	300 µl of B7-H3/CD276 (0.625 ng /ml)	300 µl of ELISA Buffer 1X
0 ng/ml	300 µl of ELISA Buffer 1X	Empty tube

8.2. Sample collection, storage and dilution

Serum : Use a serum separator tube. Let samples clot at room temperature for 30 minutes before centrifugation for 20 minutes at 1,000xg. Assay freshly prepared serum or store serum in aliquot at ≤ -20°C for later use. Avoid repeated freeze/thaw cycles.

Plasma : Collect plasma using heparin, citrate or EDTA as an anticoagulant. Centrifuge for 15 minutes at 1000xg within 30 minutes of collection. Assay freshly prepared plasma or store plasma sample in aliquot at ≤ -80°C for later use. Avoid repeated freeze/ thaw cycles.

Serum, Plasma and Cell Culture Supernatant have to be diluted in ELISA Buffer 1X. Samples containing visible precipitates must be clarified before use.

NOTE: As a starting point, 1/200 dilution of serum or of plasma is recommended! If sample values fall outside the detection range of the assay, a lower or higher dilution may be required!

8.3. Assay Procedure (Checklist)

☐	1. Determine the number of 16-well strips needed for the assay and insert them in the frame for current use. The extra strips are left in the bag with 2 silica gel minibags and stored at 4°C. NOTE: <i>Remaining 16-well strips coated with B7-H3/CD276 antibody when opened can be stored in the presence of 2 silica gel minibags at 4°C for up to 1 month.</i>
☐	2. Add 100 µl of the different standards into the appropriate wells in duplicate! At the same time, add 100 µl of diluted plasma, serum or cell culture supernatant samples in duplicate to the wells (see 8.1. Preparation and Storage of Reagents and 8.2 Preparation of Samples).
☐	3. Cover the plate with plastic film and incubate for 2 hours at Room Temperature.
☐	4. Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of five washes. After the last wash, complete removal of liquid is essential for good performance.
☐	5. Add 100 µl to each well of the diluted Detection Antibody (DET) (see 8.1 Preparation and Storage of Reagents).
☐	6. Cover the plate with plastic film and incubate for 1 hour at Room Temperature.
☐	7. Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of five washes. After the last wash, complete removal of liquid is essential for good performance.
☐	8. Add 100 µl to each well of the diluted HRP Labeled Streptavidin (STREP-HRP) (see 8.1. Preparation and Storage of Reagents).
☐	9. Cover the plate with plastic film and incubate for 30 min at Room Temperature.
☐	10. Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of five washes. After the last wash, complete removal of liquid is essential for good performance.
☐	11. Add 100 µl to each well of TMB substrate solution (TMB).
☐	12. Allow the color reaction to develop at Room Temperature in the dark for 10-15 minutes. Do not cover the plate.
☐	13. Stop the reaction by adding 100 µl of Stop Solution (STOP). Tap the plate gently to ensure thorough mixing. The substrate reaction yields a blue solution that turns yellow when Stop Solution (STOP) is added.
	! CAUTION: CORROSIVE SOLUTION !
☐	14. Measure the OD at 450 nm in an ELISA reader.

9. Calculation of Results

- Average the duplicate readings for each standard and sample and subtract the average blank value (obtained with the 0 ng/ml point).
- Generate the standard curve by plotting the average absorbance obtained for each standard concentration on the horizontal (X) axis vs. the corresponding B7-H3/CD276 concentration (ng/ml) on the vertical axis (see **10. TYPICAL DATA**).
- Calculate the B7-H3/CD276 concentrations of samples by interpolation of the regression curve formula as shown above in a form of a quadratic equation
- If the test sample was diluted, multiply the interpolated value by the dilution factor to calculate the concentration of human B7-H3/CD276 in the sample.

10. Typical Data

The following data are obtained using the different concentrations of standard as described in this protocol:

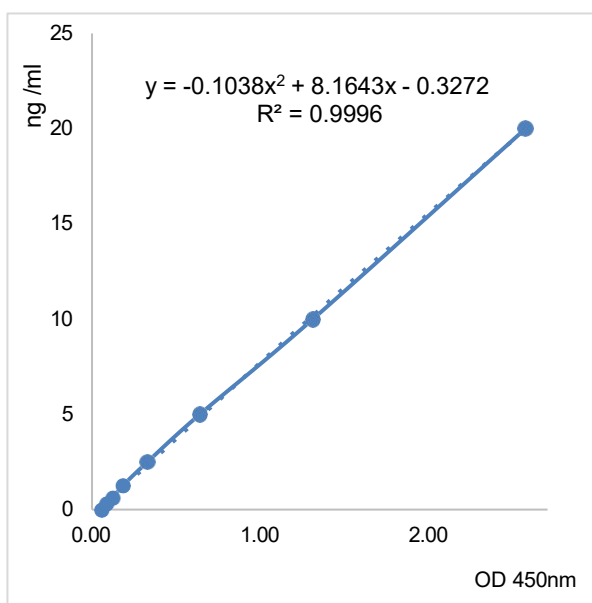


Figure: Standard curve

Standard B7-H3/CD276 (ng/ml)	Optical Density (mean)
20	2.570
10	1.309
5	0.639
2.5	0.329
1.25	0.182
0.625	0.123
0.3125	0.087
0	0.058

11. Performance Characteristics

A. Sensitivity (Limit of detection):

The lowest level of human B7-H3/CD276 that can be detected by this assay is **<0.3 ng/ml**.

NOTE: The Limit of detection was measured by adding three standard deviations to the mean value of 30 zero standard replicates.

B. Assay range: 0.3125 ng/ml – 20 ng/ml

C. Specificity:

This ELISA is specific for the measurement of natural and recombinant human B7-H3/CD276.

D. Intra-assay precision:

Four samples of known concentrations of human B7-H3/CD276 were assayed in replicates 4 times to test precision within an assay.

Samples	Means (ng/ml)	SD	CV (%)	n
A1	951.18	10.64	1.12	4
A2	697.51	54.96	7.88	4
A3	782.57	29.85	3.81	4
A4	1172.33	74.49	6.35	4

E. Inter-assay precision:

Four samples of known concentrations of human B7-H3/CD276 were assayed in 4 separate assays to test precision between assays.

Samples	Means (ng/ml)	SD	CV (%)	n
B1	1071.76	33.12	3.09	4
B2	1046.14	65.62	6.27	4
B3	733.66	26.83	3.66	4
B4	1593.98	91.76	5.76	4

F. Recovery:

When samples are spiked with known concentrations of human B7-H3/CD276, the recovery averages range from 85% to 97%.

G. Linearity:

Different samples containing human B7-H3/CD276 were diluted several folds (1/200 to 1/800 for sera and plasmas) and the measured recoveries ranged from 88% to 118%.

H. Expected values:

Human B7-H3/CD276 levels range in serum or plasma from <500 ng/ml to >2'000 ng/ml.

12. Technical Hints and Limitations

- It is recommended that all standards and samples be run in duplicate.
- Do not combine leftover reagents with those reserved for additional wells.
- Reagents from the kit with a volume less than 100µl should be centrifuged.
- Residual wash liquid should be drained from the wells after last wash by tapping the plate on absorbent paper.
- Crystals could appear in the 10X solution due to high salt concentration in the stock solutions. Crystals are readily dissolved at room temperature or at 37°C before dilution of the buffer solutions.
- Once reagents have been added to the 16-well strips, DO NOT let the strips DRY at any time during the assay.
- Keep TMB Solution protected from light.
- The Stop Solution (STOP) consists of sulfuric acid. Although diluted, the Stop Solution should be handled with gloves, eye protection and protective clothing.

13. Troubleshooting

PROBLEM	POSSIBLE CAUSES	SOLUTIONS
No signal or weak signal	Omission of key reagent	Check that all reagents have been added in the correct order.
	Washes too stringent	Use an automated plate washer if possible.
	Incubation times inadequate	Incubation times should be followed as indicated in the manual.
	Plate reader settings not optimal	Verify the wavelength and filter setting in the plate reader.
	Incorrect assay temperature	Use recommended incubation temperature. Bring substrates to room temperature before use.
High background	Concentration of STREP-HRP too high	Use recommended dilution factor.
	Inadequate washing	Ensure all wells are filling wash buffer and are aspirated completely.
Poor standard curve	Wells not completely aspirated	Completely aspirate wells between steps.
	Reagents poorly mixed	Be sure that reagents are thoroughly mixed.
Unexpected results	Omission of reagents	Be sure that reagents were prepared correctly and added in the correct order.
	Dilution error	Check pipetting technique and double-check calculations.

14. Notes