

## PCSK9(D374T) [Biotinylated]-LDLR Binding Assay Kit

**Description**

The PCSK9(D374T) [Biotinylated]-LDLR Binding Assay Kit is designed for screening and profiling purposes. The kit comes in a convenient 384-well format, with biotin-labeled mutant protein PCSK9(D374T), purified LDLR ectodomain, streptavidin-labeled HRP, and assay buffer for 384 reactions. This assay takes advantage of the high sensitivity of detection of biotin-labeled PCSK9(D374T) by streptavidin-HRP. Only a few steps are required for the assay. First, LDLR ectodomain is coated on a 384-well plate. Next, PCSK9(D374T) is incubated with LDLR on the plate. Finally, the plate is treated with streptavidin-HRP followed by addition of an HRP substrate to produce chemiluminescence, which can then be measured using a chemiluminescence reader.

**Background**

PCSK9 (Proprotein convertase subtilisin/kexin type 9) functions as a negative regulator of hepatic low-density lipoprotein receptors (LDLRs) by binding to the LDLR ectodomain. The D374T mutation is associated with hypercholesterolemia; this form of PCSK9 is more potent at decreasing LDL uptake than wild-type PCSK9, most likely by increasing the binding affinity of PCSK9 for the LDLR.

**Applications**

Study enzyme kinetics and screen small molecular inhibitors in high throughput (HTS) applications.

**Supplied Materials**

Catalog #	Name	Amount	Storage	
71211	PCSK9(D374T), Biotin-Labeled*	2 x 10 µg	-80°C	<b>Avoid multiple freeze/thaw cycles</b>
71205	LDLR, FLAG-Tag*	2 x 10 µg	-80°C	
79742	Streptavidin-HRP	2 x 10 µl	4°C	
33298	3x PL-01 Assay Buffer	2 x 50 ml	-20°C	
79728	Blocking Buffer 2	2 x 50 ml	4°C	
79670	ELISA ECL Substrate A (translucent bottle)	2 x 6 ml	Room Temp	
	ELISA ECL Substrate B (brown bottle)	2 x 6 ml	Room Temp	
78188	384-well plate	1	Room Temp	

\*The concentration of the protein is lot-specific and will be indicated on the tube.

**Materials Required but Not Supplied**

- PBS buffer (Phosphate Buffer Saline)
- Luminometer or microplate reader capable of reading chemiluminescence
- Rotating or rocker platform

**Storage Conditions**

This assay kit will perform optimally for up to 6 months from date of receipt when the materials are stored as directed. **Avoid multiple freeze/thaw cycles!**

**Safety**

This product is for research purposes only and not for human or therapeutic use. This product should be considered hazardous and is harmful by inhalation, in contact with skin, eyes, clothing, and if swallowed. If contact occurs, wash thoroughly.

**Assay Protocol**

- All samples and controls should be performed in duplicates
- The assay should include a “Blank”, a “Positive control” and a “Negative control”

**Coating the plate with LDLR:**

- 1) Thaw LDLR on ice. Briefly spin the tube containing LDLR to recover the full contents of the tube. Aliquot into single use aliquots depending on how many times the plate will be used and immediately store the aliquots at -80°C.
- 2) Dilute LDLR to 2 ng/μl (25 μl/well) in PBS.
- 3) Add 25 μl of diluted LDLR solution to each well and incubate overnight at 4°C. Leave a couple of wells empty (uncoated), for use as a “Negative Control”.
- 4) The following day, dilute 3x PL-01 Assay Buffer to 1x PL-01 Assay Buffer by adding one volume of 3x PL-01 Assay Buffer to two volumes of distilled water.
- 5) Remove LDLR solution from the plate. Wash the plate 3 times with 50 μl/well of 1x PL-01 Assay Buffer. Tap the plate onto clean paper towels to remove the liquid.
- 6) Block by adding 50 μl of Blocking Buffer 2 to each well. Incubate for 1 hour at room temperature.
- 7) Remove the Blocking Buffer.

**Step 1:**

- 1) Add 12.5 μl of 1x PL-01 Assay buffer to the “Positive control”, “Test Inhibitor”, and “Negative Control”. Add 22.5 μl of 1x PL-01 Assay buffer to the “Blank”
- 2) Thaw PCSK9(D374T) on ice. Briefly spin the tube containing the enzyme to recover the full contents of the tube. Aliquot PCSK9(D374T) into single use aliquots depending on how many times the plate will be used and immediately store the aliquots at -80°C
- 3) Dilute PCSK9(D374T) in 1x PL-01 Assay Buffer at 2.5 ng/μl (25 ng/10 μl/well). Keep the diluted protein on ice until use. Discard any unused diluted protein after use.

	<b>Blank</b>	<b>Negative Control</b>	<b>Positive Control</b>	<b>Test Inhibitor</b>
1x PL-01 Assay Buffer	22.5 μl	12.5 μl	12.5 μl	12.5 μl
Diluent Solution	2.5 μl	2.5 μl	2.5 μl	–
Test inhibitor/activator	–	–	–	2.5 μl
PCSK9(D374T) (2.5 ng/μl)	–	10 μl	10 μl	10 μl
<b>Total</b>	<b>25 μl</b>	<b>25 μl</b>	<b>25 μl</b>	<b>25 μl</b>

- 4) Prepare the compound.
  - a. If the compound is dissolved in DMSO, prepare a solution at a concentration 100-fold higher than the final desired concentration. Then dilute 10-fold in 1x PL-01 Assay Buffer.

*Note: To run an  $IC_{50}$  or test various concentrations of the compound, prepare serial dilutions using PL-01 Assay Buffer containing 10% DMSO, so the final concentration of DMSO will be 1% in all samples.*

- b. If the compound is soluble in water, prepare a solution of the compound in 1x PL-01 Assay Buffer that is 10-fold higher than the final desired concentration.
- 5) Add 2.5  $\mu$ l of inhibitor solution to each well designated "Test Inhibitor." For the "Positive Control," "Negative Control" and "Blank," add 2.5  $\mu$ l of diluent solution (for example, 10% DMSO in water if the inhibitor was dissolved in DMSO).
- 6) Initiate the reaction by adding 10  $\mu$ l of diluted PCSK9(D374T) to all wells except the "Blank".
- 7) Incubate at room temperature for two hours.
- 8) Remove the reaction solution and wash the plate 3 times with 50  $\mu$ l/well of 1x PL-01 Assay Buffer. Tap the plate onto clean paper towels to remove the liquid.
- 9) Block by adding 50  $\mu$ l of Blocking Buffer 2 to each well. Incubate for 10 minutes at room temperature. Decant to remove supernatant.

#### **Step 2:**

- 1) Dilute Streptavidin-HRP 1000-fold with Blocking Buffer 2.
- 2) Add 50  $\mu$ l to each well. Incubate for 1 hour at room temperature with slow shaking.
- 3) Wash the plate three times with 50  $\mu$ l/well of 1x PL-01 Assay Buffer. Tap the plate onto clean paper towels to remove the liquid.
- 4) Block by adding 50  $\mu$ l of Blocking Buffer 2 to each well. Incubate for 10 minutes at room temperature. Remove the Blocking Buffer and tap the plate onto clean paper towels to remove the liquid.
- 5) Just before use, mix 25  $\mu$ l/well of ELISA ECL substrate A and 25  $\mu$ l/well of ELISA ECL substrate B, then add 50  $\mu$ l to each well. Discard any unused chemiluminescent reagent after use.
- 6) Immediately read the plate in a luminometer or plate reader capable of reading chemiluminescence. The "Blank" value is subtracted from all readings.

#### **Reading Chemiluminescence**

Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry. To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second, delay after plate movement is 100 msec. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the "hole" position on the filter wheel; Optics position: Top; Read type: endpoint. Sensitivity may be adjusted based on the luminescence of a control assay without enzyme (Reader was set at 100 for validation data).

## Example Results

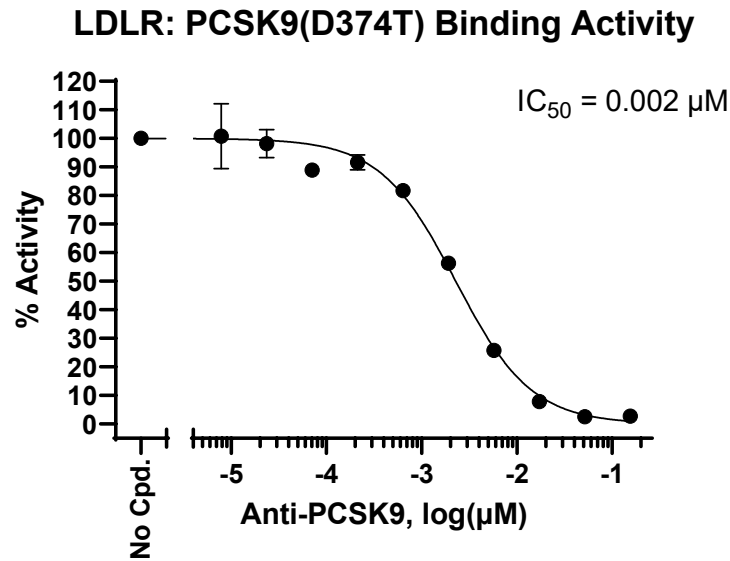


Figure 1. PCSK9-LDLR binding activity.

Binding of PCSK8 to LDLR was measured using the PCSK9(D374T) [Biotinylated]-LDLR Binding Assay Kit (BPS Bioscience #78326) with increasing concentrations of Anti-PCSK9 Neutralizing Antibody (BPS Bioscience #71207). Luminescence was measured using a Bio-Tek fluorescent microplate reader.

Data shown is representative. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com).

### Troubleshooting Guide

Visit [bpsbioscience.com/assay-kits-faq](https://bpsbioscience.com/assay-kits-faq) for detailed troubleshooting instructions. For all further questions, please email [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

### Related Products:

Products	Catalog #	Size
PCSK9, His-tag Recombinant	71204	50 $\mu\text{g}$
PCSK9, C-terminal His-Avi-Tag, Biotin-Labeled Recombinant	71304	20 $\mu\text{g}$
LDLR, Biotin-labeled Recombinant	71206	20 $\mu\text{g}$
Anti-PCSK9 Neutralizing Antibody	71207	50 $\mu\text{g}$
PCSK9-LDLR TR-FRET Assay Kit	72010	384 rxns.
PCSK9(D374T)-LDLR TR-FRET Assay Kit	72011	384 rxns.