

## Product datasheet

# Paraoxonase 1 Activity Assay Kit ab241044

[3 Images](#)

### Overview

<b>Product name</b>	Paraoxonase 1 Activity Assay Kit
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Serum, Plasma
<b>Product overview</b>	Paraoxonase 1 Activity Assay Kit (ab241044) enables rapid measurement of Paraoxonase 1 (PON1) activity, utilizing a fluorogenic substrate that is converted into a highly fluorescent product (Ex/Em = 368/460 nm). This ensures dramatically greater sensitivity than UV or colorimetric assays and eliminates the need for dangerous toxic substrates. A selective PON1 inhibitor is provided for verification of PON1 specific activity. The assay is simple to perform, high-throughput adaptable and can detect a minimum of 2.0 $\mu$ U paraoxonase activity with a sample volume of 5 $\mu$ L.
<b>Platform</b>	Microplate reader

### Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	100 tests
Fluorescence Standard	1 vial
Paraoxonase Assay Buffer	1 x 50ml
Paraoxonase Positive Control	1 vial
PON1 Inhibitor (2-hydroxyquinoline)	1 vial
PON1 Substrate	1 vial

**Function** Hydrolyzes the toxic metabolites of a variety of organophosphorus insecticides. Capable of hydrolyzing a broad spectrum of organophosphate substrates and lactones, and a number of aromatic carboxylic acid esters. Mediates an enzymatic protection of low density lipoproteins against oxidative modification and the consequent series of events leading to atheroma formation.

**Tissue specificity** Plasma, associated with HDL (at protein level). Expressed in liver, but not in heart, brain,

placenta, lung, skeletal muscle, kidney or pancreas.

### Involvement in disease

Genetic variation in PON1 is associated with susceptibility to microvascular complications of diabetes type 5 (MVCD5) [MIM:612633]. These are pathological conditions that develop in numerous tissues and organs as a consequence of diabetes mellitus. They include diabetic retinopathy, diabetic nephropathy leading to end-stage renal disease, and diabetic neuropathy. Diabetic retinopathy remains the major cause of new-onset blindness among diabetic adults. It is characterized by vascular permeability and increased tissue ischemia and angiogenesis. Note=Homozygosity for the Leu-54 allele is strongly associated with the development of retinal disease in diabetic patients.

### Sequence similarities

Belongs to the paraoxonase family.

### Post-translational modifications

Glycosylated.

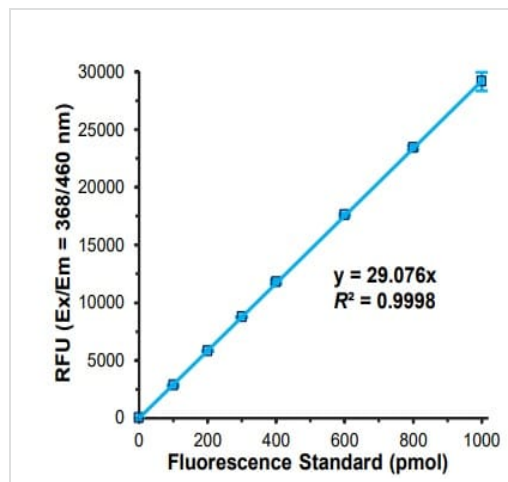
The signal sequence is not cleaved.

Present in two forms, form B contains a disulfide bond, form A does not.

### Cellular localization

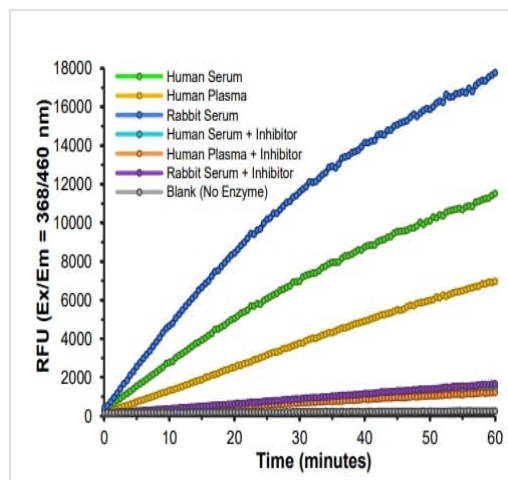
Secreted > extracellular space.

## Images



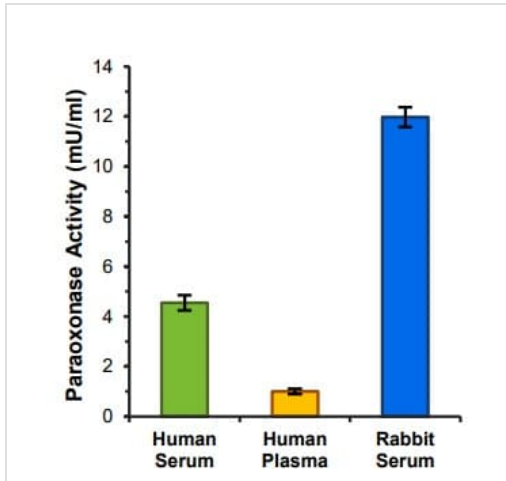
Standard curve of PON1 Substrate metabolite fluorescence. One mole of fluorescence standard corresponds to the metabolism of one mole of PON1 Substrate.

Standard curve of PON1.



Reaction kinetics of PON1 Substrate metabolism in donor-pooled human serum (5  $\mu$ L), donor-pooled human plasma (5  $\mu$ L) and rabbit serum (2.5  $\mu$ L) in the presence and absence of 200  $\mu$ M of the selective PON1 inhibitor 2- hydroxyquinoline (no inhibitor conditions contained 0.4% DMSO as a solvent control).

Reaction kinetics of PON1 Substrate metabolism.



Quantification of PON1 activity in serum/plasma samples (mean  $\pm$  SEM of four independent replicates). Assays were performed according to the kit protocol.

Quantification of PON1 activity in serum/plasma samples.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors