

C-Peptide ELISA Kit ab178641

3 References 1 Image

Overview

Product name	C-Peptide ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	C-peptide	16		
				CV%
				<=6.2%
	Inter-assay			
	Sample	n	Mean	SD
	C-peptide	20		
				CV%
				<=10%
Sample type	Serum, Hep Plasma, Cit plasma			
Assay type	Sandwich (quantitative)			
Sensitivity	0.01 ng/ml			
Range	0.7 ng/ml - 1.9 ng/ml			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	C-peptide ELISA kit (ab178641) is designed for the accurate quantitative measurement of C-peptide in Human serum and plasma.			

A 96-well plate has been precoated with Streptavidin. Samples, standards and the C-peptide HRP and Biotin conjugate are added to the wells. Biotinylated monoclonal and horseradish peroxidase (HRP) labelled antibodies are added and the reactants are mixed. The different types of antibodies used have high affinity and specificity and are directed against distinct and different epitopes of C-Peptide. Reaction between the various C-Peptide antibodies and native C-Peptide occurs in the microwells without competition or steric hindrance forming a soluble sandwich complex. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is directly proportional to the amount of C-peptide in the sample and the intensity is measured at 450 nm.

Notes	C-peptide is the abbreviation for connecting peptide; it is a 31-aminoacid peptide. C-peptide of insulin is the C-terminal cleavage product produced during processing of the insulin pro-hormone to the mature insulin molecule. Proinsulin is cleaved when it is released from the pancreas into the blood - one C-peptide for each insulin molecule. C-Peptide is devoid of any biological activity but appears to be necessary to maintain the structural integrity of Insulin.
Platform	Microplate

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
50X Washing Solution	1 x 20ml
Cover Foil	1 unit
C-Peptide HRP and Biotin Conjugate	1 x 13ml
C-Peptide Standard 0 - 0 ng/mL (Lyophilized)	1 vial
C-Peptide Standard 1 - 0.2 ng/mL (Lyophilized)	1 vial
C-Peptide Standard 2 - 1.0 ng/mL (Lyophilized)	1 vial
C-Peptide Standard 3 - 2.0 ng/mL (Lyophilized)	1 vial
C-Peptide Standard 4 - 5.0 ng/mL (Lyophilized)	1 vial
C-Peptide Standard 5 - 10.0 ng/mL (Lyophilized)	1 vial
Stop Solution	1 x 15ml
Streptavidin Coated Microplate (12 x 8 wells)	1 unit
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

Function Insulin decreases blood glucose concentration. It increases cell permeability to monosaccharides, amino acids and fatty acids. It accelerates glycolysis, the pentose phosphate cycle, and glycogen synthesis in liver.

Involvement in disease Defects in INS are the cause of familial hyperproinsulinemia (FHPRI) [MIM:176730]. Defects in INS are a cause of diabetes mellitus insulin-dependent type 2 (IDDM2) [MIM:125852]. IDDM2 is a multifactorial disorder of glucose homeostasis that is characterized by susceptibility to ketoacidosis in the absence of insulin therapy. Clinical features are polydipsia, polyphagia and polyuria which result from hyperglycemia-induced osmotic diuresis and secondary thirst. These derangements result in long-term complications that affect the eyes, kidneys, nerves, and blood vessels. Defects in INS are a cause of diabetes mellitus permanent neonatal (PNDM) [MIM:606176]. PNDM is a rare form of diabetes distinct from childhood-onset autoimmune diabetes mellitus type

1. It is characterized by insulin-requiring hyperglycemia that is diagnosed within the first months of life. Permanent neonatal diabetes requires lifelong therapy.

Defects in INS are a cause of maturity-onset diabetes of the young type 10 (MODY10) [MIM:613370]. MODY10 is a form of diabetes that is characterized by an autosomal dominant mode of inheritance, onset in childhood or early adulthood (usually before 25 years of age), a primary defect in insulin secretion and frequent insulin-independence at the beginning of the disease.

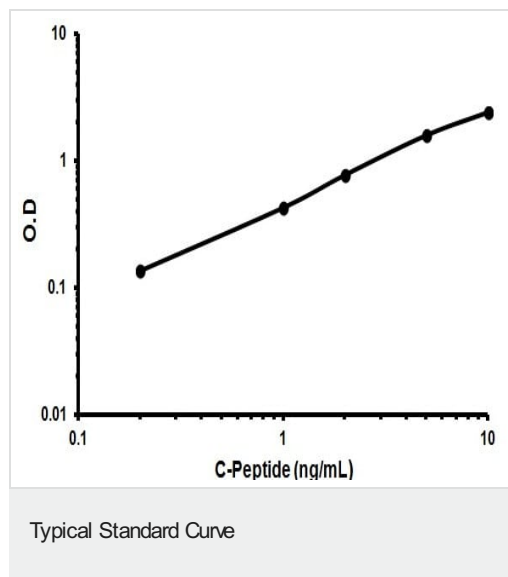
Sequence similarities

Belongs to the insulin family.

Cellular localization

Secreted.

Images



Representative standard curve using ab178641.

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