

# Product Information

sigma-aldrich.com

3050 Spruce Street, Saint Louis, MO 63103 USA  
Tel: (800) 521-8956 (314) 771-5765 Fax: (800) 325-5052 (314) 771-5757  
email: techservice@sial.com sigma-aldrich.com

## Carboxypeptidase A Assay Kit

Catalog Number **CS1130**

Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Carboxypeptidase A (CPA) catalyzes the hydrolysis of C-terminal aromatic or aliphatic amino acids of proteins or peptides. Carboxypeptidase A is found in the pancreas and in mast cells. Carboxypeptidase A found in mast cells (MC-CPA) is a highly conserved secretory granule protease that is important for mast cell differentiation.<sup>1</sup> MC-CPA levels rise during anaphylaxis, a severe acute allergic reaction.<sup>2</sup> This enzyme is considered to be a sensitive marker for anaphylaxis, much more sensitive than the present marker tryptase. It has been suggested that MC-CPA can degrade snake venom components and honeybee venoms, thus, having a protective function during envenomation.<sup>3</sup>

The kit allows the fast and convenient determination of CPA activity in biological samples, as well as screening for enzyme inhibitors. The kit contains all the necessary components including a Carboxypeptidase A control enzyme and Carboxypeptidase Inhibitor.

The CPA activity assay is based on the hydrolysis of the substrate N-(4-methoxyphenylazoformyl)-Phe-OH. The substrate has a strong absorption peak at 350 nm ( $\epsilon^{\text{mM}} = 19.0$ ). The reaction progression is reflected by a decrease in absorption at 350 nm.<sup>4</sup>

Unit definition: 1 unit will hydrolyze 1  $\mu\text{mole}$  of N-(4-methoxyphenylazoformyl)-Phe-OH per minute at pH 8 at  $25\text{ }^{\circ}\text{C}$ .

### Components

The kit is sufficient for 200 assays in 96 well plates or 40 assays in 1 mL cuvette.

Assay Buffer 5 mL  
Catalog Number A7731

Substrate Solution 50  $\mu\text{L}$   
Catalog Number S8697

Carboxypeptidase A 200  $\mu\text{L}$   
from bovine pancreas  
Catalog Number C8368

Carboxypeptidase Inhibitor 200  $\mu\text{L}$   
from potato tuber

Catalog Number C8493

Sodium Carbonate 2.5 g  
Catalog Number S2127

### Reagents and Equipment Required but Not Provided

- 96 well clear flat bottom plate (Catalog Number CLS3358)
- Plate reader or spectrophotometer
- 1 mL quartz cuvette (optional)
- Ultrapure water
- Magnetic stirrer and stir bar

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Preparation Instructions

Use ultrapure water (17 M $\Omega$ -cm or equivalent) for the preparation of reagents and throughout the procedure.

Stop Solution - Add 24 mL of ultrapure water to the sodium carbonate bottle (Catalog Number S2127) and mix well on a magnetic stirrer until completely dissolved. Store the Stop Solution at room temperature.

Reaction Mixture – Prepare 1 mL of reaction mixture by mixing 800  $\mu\text{L}$  of ultrapure water, 200  $\mu\text{L}$  of Assay Buffer (Catalog Number A7731), and 2  $\mu\text{L}$  of the Substrate Solution (Catalog Number S8697). One mL of Reaction Mixture is sufficient for 2 assays in 1 mL cuvette or 10 assays in a 96 well plate.

### Storage/Stability

The kit is shipped on dry ice and storage at  $-20\text{ }^{\circ}\text{C}$  is recommended. Upon arrival store the Sodium Carbonate (Catalog Number S2127) at room temperature.

### Procedure

The following procedure is for 96 well plate assays. Perform the reactions in duplicates.

**Notes:** The 96 well plate assay is based on a fast end-point reaction. Therefore, if several reactions are performed simultaneously, it is highly recommended to use a multichannel pipette for initiation and termination of the reaction.

It is possible to perform the assay in cuvettes. The volumes of the reagents should be adjusted according to the increase in the reaction volume.

**Table 1.**

Reaction scheme for 96 well plate assay

Sample	Enzyme	Inhibitor	Water	Reaction Mixture
Blank	–	–	100 $\mu$ L	100 $\mu$ L
Test sample	x $\mu$ L sample	–	100–x $\mu$ L	100 $\mu$ L
Positive Control	1–2 $\mu$ L Carboxypeptidase A	–	97–98 $\mu$ L	100 $\mu$ L
Control with Inhibitor	2 $\mu$ L Carboxypeptidase A	1 $\mu$ L	97 $\mu$ L	100 $\mu$ L

x - volume of the test sample added to the reaction

1. Prepare a blank control by adding 100  $\mu$ L of ultrapure water to the appropriate wells.
2. Prepare a test sample reaction by adding 1–50  $\mu$ L (1–100  $\mu$ g protein) of sample to the appropriate wells. Adjust the volume to 100  $\mu$ L with ultrapure water.
3. Prepare a positive control by adding 1–2  $\mu$ L of the Carboxypeptidase A control enzyme to the appropriate wells. Adjust the volume to 100  $\mu$ L with ultrapure water.
4. Prepare a control with inhibitor reaction by adding 2  $\mu$ L of the Carboxypeptidase A control enzyme and 1  $\mu$ L of Carboxypeptidase Inhibitor to the appropriate wells. Adjust the total volume to 100  $\mu$ L with ultrapure water.

**Note:** If other enzyme preparations or/and inhibitors are used, their volumes must be optimized based on their activities. Add to a well the appropriate amounts of inhibitor and enzyme of choice, and adjust the total volume to 100  $\mu$ L with ultrapure water. Then continue with the procedure.

5. Start the reaction by the addition of 100  $\mu$ L of Reaction Mixture to each well.
6. Incubate the plate for 3–5 minutes at 25  $^{\circ}$ C.
7. Stop the reaction with 100  $\mu$ L of Stop Solution.  
**Note:** Some detergents (e.g., TRITON<sup>®</sup> X-100) may precipitate when the sodium carbonate Stop Solution is added. If a precipitate forms, measure the absorption without the addition of Stop Solution.
8. Read the absorption at 350 nm.

### Results

#### Calculation

The reaction rate is linear only when the decrease in absorption ( $A_{350}$ ) does not exceed 0.35. For calculating the CPA activity according to the following equations, use only the linear range of the reaction.

Calculation of the CPA activity detected using 96 well plate:

$$\text{CPA Activity (milliunits/mL)} = \frac{\Delta A_{350} \times \text{Dilution} \times 0.2}{\text{Time} \times 0.019 \times 0.55 \times V_{\text{sample}}}$$

Calculation of the CPA activity detected using a 1 mL cuvette:

$$\text{CPA Activity (milliunits/mL)} = \frac{\Delta A_{350} \times \text{Dilution}}{\text{Time} \times 0.019 \times V_{\text{sample}}}$$

Where:

$\Delta A_{350}$  - difference between initial and final absorption at 350 nm

**Note:** initial absorption is equal to the blank control absorption

**Dilution** - dilution of the test sample or enzyme stock

**Time** - reaction duration (incubation time) from the beginning of the reaction to the end (minutes)

$V_{\text{sample}}$  - volume of the sample/control enzyme added to the reaction mixture (mL)

**0.019** -  $\mu$ molar extinction coefficient of the substrate

**0.2** - volume of reaction in mL

**0.55** - effective pathlength in 96 well plates

### Calibration curve

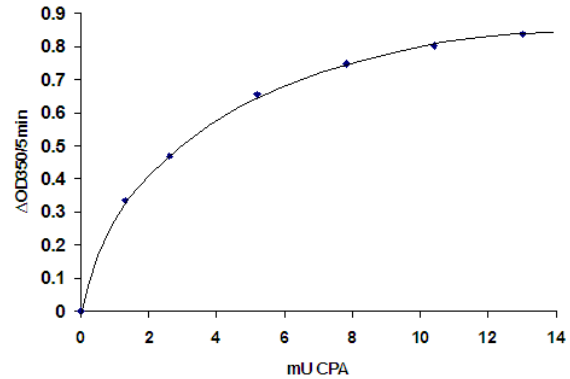
The calibration curve for the supplied Carboxypeptidase A control is not linear (see Figure 1). Therefore, to determine the activity of an unknown sample, which is not in the linear range of the reaction, a calibration curve should be used. The activity of the unknown sample may be determined by this curve.

Determination of a Carboxypeptidase A calibration curve:

1. Dilute 40  $\mu\text{L}$  of the Carboxypeptidase A control enzyme (Catalog Number C8368) 10-fold with ultrapure water.
2. Place 0, 10, 20, 40, 60, and 80  $\mu\text{L}$  of the diluted Carboxypeptidase A control enzyme into the appropriate wells. Bring the volume to 100  $\mu\text{L}$  with ultrapure water.
3. Start the reaction by the addition of 100  $\mu\text{L}$  of Reaction Mixture.
4. Incubate the plate for 5 minutes at 25  $^{\circ}\text{C}$ .
5. Stop the reaction with 100  $\mu\text{L}$  of Stop Solution.
6. Read the absorbance at 350 nm.
7. Plot the calibration curve ( $\Delta\text{OD}_{350}$  per minute as a function of CPA activity, see Figure 1).

**Note:** CPA activity of the control enzyme must be calculated according to the equation in the Results section, using a single point where the reaction rate is still in the linear range: Use the  $\Delta\text{OD}_{350}$  ( $\Delta A_{350}$ ) value for the first point (10  $\mu\text{L}$ ) of the control enzyme in the equation and determine the activity in milliunits. Calculate the activity for the rest of the enzyme calibration curve samples according to the value obtained for the 10  $\mu\text{L}$  sample (e.g., if there are 0.9 milliunits in the 10  $\mu\text{L}$  sample, then the 20  $\mu\text{L}$  sample contains 1.8 milliunits).

**Figure 1.**  
Carboxypeptidase A Calibration Curve



Calibration Curve for the control Carboxypeptidase A provided with this kit. It is possible that a slightly different curve may be obtained for CPA from other biological sources.

### References

1. Feyerabend, T.B., et al., Loss of histochemical identity in mast cells lacking Carboxypeptidase A. *Mol. Cell Biol.*, **25**, 6199–6210 (2005).
2. Zhou, X., et al., Mast cell Carboxypeptidase as a new clinical marker for anaphylaxis. *J. Allergy Clin. Immunol.*, **117**, S85 (2006).
3. Metz, M., et al., Mast cells can enhance resistance to snake and honeybee venoms. *Science*, **313**, 526-530 (2006).
4. Mock, W.L., et al., Arazoformyl peptide surrogates as spectrophotometric kinetic assay substrates for Carboxypeptidase A. *Anal. Biochem.*, **239**, 218-222 (1996).

TRITON is a registered trademark of Union Carbide Corp.

EB, MAM 01/08-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.