



Caspase Substrates & Set

INTRODUCTION

Caspase Substrates & Sets are ready-to-use solutions of caspase substrates for both fluorometric and colorimetric detection of caspase activity in various cell types. These substrates are prepared as 1mM solution and are supplied individually as well as in a set of all caspase substrates for studying multiple apoptosis processes.

ITEM(S) SUPPLIED

Catalog #	Substrate	Size
CPS-145	Caspase-1,4,5 Substrate; Ac-WEHD-AFC (1mM)	250µl
CPS-002	Caspase-2 Substrate; Ac-VDVAD-AFC (1mM)	250µl
CPS-370	Caspase-3,7,10 Substrate; Ac-DEVD-AFC (1mM)	250µl
CPS-006	Caspase-6 Substrate; Ac-VEID-AFC (1mM)	250µl
CPS-008	Caspase-8 Substrate; Ac-IETD-AFC (1mM)	250µl
CPS-009	Caspase-9 Substrate; Ac-LEHD-AFC (1mM)	250µl
CPS-013	Caspase-13 Substrate; Ac-LEED-AFC (1mM)	250µl
CPS-113S	Caspase 1-10 and 13 <u>Substrates Set</u>	250µl each

STORAGE CONDITION

It is shipped in blue ice. Store at -20°C and is stable for 6-8 months if stored properly.

PROTOCOL

PREPARATION BEFORE USE

Take out the vial(s) from refrigerator and thaw to RT. Vortex to mix the contents. Avoid repeated freezing and thawing (prepare small aliquots as per requirement). The substrate is ready to use for apoptosis screening and other related assays.

RELATED PRODUCTS:

CasPASE™-Apoptosis Assay Kit (Cat # 786-200 to 786-206)

The assay kit provides a simple method for assaying various caspase activity. For details please visit our web site www.GBiosciences.com.

CasPASE Inhibitors and Set Ready to use caspase inhibitors are provided individually as well as in a set. For details please visit our web site www.GBiosciences.com

NOTE: For other related products, visit our web site at www.GBiosciences.com or contact us.

Rev 05.21.09/MM

