



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Fluorogenic DPP3 Assay Kit

Catalog #: 80203

DESCRIPTION: Dipeptidyl peptidase-3 (DPP3) is a serine exopeptidase that cleaves X-proline dipeptides from the N-terminus of polypeptides. DPP3 activity has been associated with endometrial and ovarian cancers. The *Fluorogenic DPP3 Assay Kit* is designed to measure DPP3 activity using purified DPP3 for screening and profiling applications. It comes in a convenient 96-well format, with purified DPP3 enzyme, DPP substrate, and DPP assay buffer for 100 enzyme reactions. The key to the *Fluorogenic DPP3 Assay Kit* is the specific, fluorogenic substrate. Using this kit, only one simple step on a microtiter plate is required for DPP3 reactions. The fluorometric substrate is incubated with a sample containing DPP3 enzyme to produce a fluorophore that can then be measured using a fluorescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
80030	DPP3 human recombinant enzyme	5 µg	-80°C	<i>Avoid freeze/thaw cycles!</i>
80300	DPP assay buffer	10 ml	-20°C	
80332	Fluorogenic DPP substrate 2 in DMSO (0.5 mM)	100 µl	-80°C	
	Fluorescent AMC standard (50 µM)	500 µl	-20°C	
	black, low binding NUNC black microtiter plate	1 plate	Room temp.	

APPLICATIONS: Great for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

REFERENCES:

1. Gevaert, K. et al., *Proteomics* **5 (14)**: 3589-3599 (2005).
2. Abramic, M. et al., *Int. J. Biochem. Cell Biol.* **36 (3)**: 434-446 (2004).

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

Immediately prior to assay:

- 1) Dilute DPP substrate 2 0.5 mM stock 5-fold with DPP buffer to make a 100 μ M solution. (Make only sufficient quantity needed for the assay; store remaining 0.5 mM stock solution in aliquots at -20°C.)
- 2) Dilute DPP3 in DPP assay buffer to 5 ng/ μ l (50 ng/reaction)*. Aliquot any remaining enzyme and store undiluted at -80°C. Keep diluted enzyme on ice. Discard any remaining diluted enzyme after use. **Note: Optimal enzyme concentration may vary with the specific activity of the enzyme.*
- 3) Dilute 25 μ l of the Fluorescent AMC standard (50 μ M stock) 2-fold with DPP buffer to make a 25 μ M solution. Make serial 2-fold dilutions of the fluorescent AMC standard in DPP buffer as follows: 12.5 μ M, 6.25 μ M, 3.12 μ M, 1.56 μ M, 0.78 μ M, 0.39 μ M, 0.19 μ M, 0.10 μ M. Aliquot the remaining 50 μ M AMC standard and store undiluted at -20°C.

Step 1:

In duplicate, add the reaction mixtures (below) to the microtiter black plate. Incubate at 22 °C for 10 min.

	Enzyme Positive Control	Test Inhibitor	AMC Standard Curve	Inhibitor Negative Control	"Blank" Negative Control
DPP3 (5 ng/ μ l)	10 μ l	10 μ l	–	–	–
DPP substrate 24 (100 μ M)	5 μ l	5 μ l	–	5 μ l	–
AMC standard (0.1 μ M – 50 μ M)	–	–	5 μ l	–	–
Inhibitor (in DPP assay buffer)	–	X μ l	–	X μ l	–
DPP assay buffer	85 μ l	85 - X μ l	95 μ l	95 - X μ l	100 μ l
Total	100 μl	100 μl	100 μl	100 μl	100 μl

Step 2:

Read sample in a microtiter-plate fluorimeter that is capable of excitation at wavelengths ranging from 350-380 nm and detection of emitted light ranging from 440-460 nm. Blank value is subtracted from all other values.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

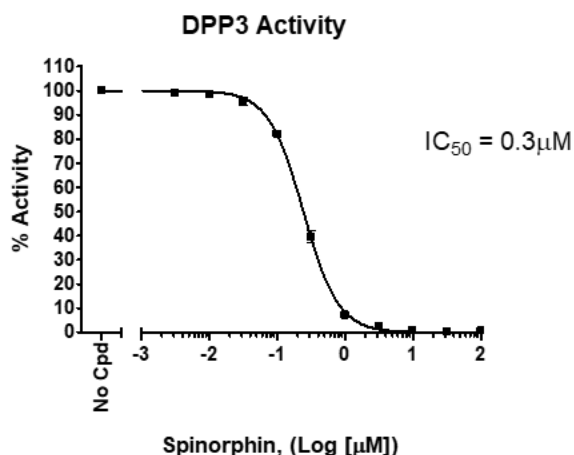
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Example of Assay Results:



DPP3 enzyme activity, measured using the *Fluorogenic DPP3 Assay Kit*, BPS Bioscience Catalog # 80203. *Note: Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com.*

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
DPP3 enzyme	80030	10 μg
DPP assay buffer	80300	20 ml
Fluorogenic DPP substrate 2	80332	100 μl
DPP4 enzyme	80040	10 μg
DPP7 enzyme	80070	10 μg
DPP8 enzyme	80080	10 μg
DPP9 enzyme	80090	10 μg
FAP enzyme	80100	10 μg
POP enzyme	80105	20 μg
Fluorogenic DPP4 assay kit	80204	96 rxns
Fluorogenic DPP7 assay kit	80207	96 rxns
Fluorogenic DPP8 assay kit	80208	96 rxns
Fluorogenic DPP9 assay kit	80209	96 rxns
Fluorogenic FAP assay kit	80210	96 rxns
Fluorogenic POP assay kit	80106	96 rxns

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Fluorogenic DPP substrate 1

80305

100 µl

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: [**info@bpsbioscience.com**](mailto:info@bpsbioscience.com)
Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)