



DPPIV Drug Discovery Kit

Designed to screen DPPIV inhibitors

Instruction Manual

BML-AK499

For research use only

✦ DPPIV Drug Discovery Kit –BML- AK499✦

BACKGROUND

DPPIV (DPP4, CD26) is a member of the class of proteases known as prolyl peptidases, which cleave proteins after proline residues¹. DPPIV, a serine dipeptidyl peptidase, cleaves the N-terminal X-Ala or X-Pro from target polypeptides, such as chemokines (e.g. CXCL11) and peptide hormones (e.g., glucagon-like peptide-1, GLP1)¹⁻³. DPPIV possesses a transmembrane region and a very short cytoplasmic domain, but is often cleaved and released as a soluble, circulating form⁴. It is found as a dimer with itself or with FAP (fibroblast activation protein- α , seprase), another prolyl peptidase¹. It also has non-peptidase functions: through its interaction with adenosine deaminase (ADA) and extracellular matrix components, it influences T-cell activation and proliferation^{2,5,6}. It is thought to play roles in diabetes, cancer, and autoimmune diseases, making it a target for drug discovery⁷⁻¹¹.

The *DPPIV Drug Discovery Kit* is a complete assay system designed to screen DPPIV inhibitors, providing enough material to perform at least 96 assays. The kit contains both a chromogenic substrate (H-Gly-Pro-pNA; $K_m=114 \mu\text{M}$) and a fluorogenic substrate (H-Gly-Pro-AMC; $K_m=50 \mu\text{M}$). Cleavage of the p-nitroaniline (pNA) from the chromogenic substrate increases absorbance at 405 nm. The fluorimetric assay is based on the cleavage of 7-amino-4-methylcoumarin (AMC) moiety from the C-terminus of the peptide substrate, which increases its fluorescence intensity at 460 nm. The kit is useful to screen inhibitors of DPPIV, a potential therapeutic target. A DPPIV inhibitor, P32/98 ($K_i=130 \text{ nM}$ ¹²), is included for use as a control.

Other DPP enzymes are available for specificity profiling. Contact Enzo Life Sciences or go to www.enzolifesciences.com.

REFERENCES:

1. J.S. Rosenblum and J.W. Kozarich *Curr. Opin. Chem. Biol.* 2003 **7** 496
2. R. Thoma *et al. Structure* 2003 **11** 947
3. H. Shibuya-Saruta *et al. J. Clin. Lab. Anal.* 1996 **10** 435
4. D.A. Pereira *et al. Braz. J. Med. Biol. Res.* 2003 **36** 567
5. I. Ben-Shooshan *et al. Biochem. Biophys. Acta* 2002 **1587** 21
6. K. Aertgeerts *et al. Protein Sci.* 2004 **13** 145
7. A. E. Weber *J. Med. Chem.* 2004 **47** 4135
8. B. Pro and N.H. Dang *Histol. Histopathol.* 2004 **19** 1345
9. U. Aytac and N.H. Dang *Curr. Drug Targets Immune Endocr. Metabol. Disord.* 2004 **4** 11
10. H. Gotoh *et al. Clin. Chem.* 1988 **34** 2499
11. D.J. Augeri *et al. J. Med. Chem.* 2005 **48** 5025
12. J.A. Pospisilik *et al. Diabetes* 2003 **52** 741

PLEASE READ ENTIRE BOOKLET BEFORE PROCEEDING WITH THE ASSAY. CAREFULLY NOTE THE HANDLING AND STORAGE CONDITIONS OF EACH KIT COMPONENT. PLEASE CONTACT ENZO LIFE SCIENCES TECHNICAL SERVICES FOR ASSISTANCE IF NECESSARY.

COMPONENTS OF BML-AK499 KIT

BML-SE434-9090 DPPIV ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant soluble human DPPIV. One U=1 μ mole/min@37°C, 100 μ M H-Gly-Pro-pNA.

STORAGE: -80°C; AVOID FREEZE/THAW CYCLES

QUANTITY: 35 mU

BML-P188-9090 pNA SUBSTRATE (H-Gly-Pro-pNA; MW=328.8)

FORM: 10 mM in DMSO

STORAGE: -80°C

QUANTITY: 150 μ L

BML-KI106-0001 pNA CALIBRATION STANDARD (p-nitroaniline; MW=138)

FORM: 50 μ M in assay buffer.

STORAGE: -80°C

QUANTITY: 1 mL

BML-P189-9090 AMC SUBSTRATE (H-Gly-Pro-AMC; MW=410.3)

FORM: 0.5 mM in DMSO

STORAGE: -80°C

QUANTITY: 150 μ L

BML-KI107-0001 AMC CALIBRATION STANDARD

(7-amino-4-methylcoumarin; MW=175)

FORM: 30 μ M in assay buffer.

STORAGE: -80°C

QUANTITY: 1 mL

BML-PI142-9090 INHIBITOR (P32/98; MW=260.4)

FORM: 1 mM in DMSO

STORAGE: -80°C

QUANTITY: 20 µL

BML-KI342-0020 ASSAY BUFFER

(50 mM Tris, pH 7.5)

FORM: Liquid in screw-cap plastic bottle

STORAGE: -80°C (or +4°C)

QUANTITY: 20 mL

80-2407 ½-VOLUME CLEAR & NBS WHITE MICROPLATE – 1 EACH

STORAGE: Ambient.

OTHER MATERIALS REQUIRED

Microplate reader capable of measuring A_{405} to ≥ 3 -decimal accuracy, or fluorescence at wavelengths of approximately 380nm (excitation)/ 460nm (emission)

Pipettes or multi-channel pipettes capable of pipetting 10-1000 µL accurately (note: dilution of reagents can be made to increase the minimal volume to >10 µL).

Ice bucket to keep reagents cold until use.

SUGGESTED EXPERIMENTAL METHODS

Note on storage: Store all components except the microplates at -80°C for the highest stability. The DPPIV enzyme should be handled carefully in order to retain maximal enzymatic activity. It is stable, in diluted or concentrated form, for several hours on ice. As supplied, DPPIV enzyme is stable for at least 5 freeze/thaw cycles. To minimize the number of freeze/thaw cycles, aliquot the DPPIV into separate tubes and store at -80°C. Do not maintain diluted components at reaction temperature (e.g. 37°C) for an extended period of time prior to running the assay.

To start assay:

1. Defrost kit components and hold on ice until use. Thaw and store DMSO components (substrates, inhibitor) at room temperature, preferably in a dark place. *Briefly centrifuge all vials.* Minimize the time that any kit component is thawed.
2. Dilute inhibitor (P32/98) 1/10 in assay buffer. Example: Add 2 μL inhibitor to 18 μL assay buffer, in a separate tube.
3. For colorimetric assay (at $A_{405\text{nm}}$, using clear microplate): Dilute substrate (H-Gly-Pro-pNA) 1/50 in assay buffer (50 μL is needed per well). Example: Add 5 μL substrate to 245 μL assay buffer, in a separate tube.
4. For fluorimetric assay (at Ex:380 nm/Em:460 nm, using white microplate), dilute substrate (H-Gly-Pro-AMC) 1/50 in assay buffer (50 μL is needed per well). Example: Add 5 μL substrate to 245 μL assay buffer, in a separate tube.
5. Add assay buffer to each desired well of microplate so that the total assay volume will be 100 μL . See **Table 1** for examples.
6. Allow microplate and diluted assay components to equilibrate to assay temperature (e.g. 37°C).
7. After a brief thawing, quickly centrifuge the vial of DPPIV (BML-SE434-9090) to bring contents to bottom of tube. Dilute enough DPPIV in assay buffer to a stock concentration of 17.3 $\mu\text{U}/\mu\text{L}$ to produce required quantity for the experiment.
8. Add DPPIV to the “Control”, “Inhibitor” and “Test Sample” wells, please note that the final amount of DPPIV *per well* is to be 0.26mU. DO NOT ADD DPPIV TO BLANKS!
9. Add 10 μL P32/98 inhibitor (diluted in step 2) to the “Inhibitor” well only! Final inhibitor concentration is 10 μM .
10. Add desired volume of test sample(s) to appropriate well(s). See **Table 1**.
11. Incubate plate for 10 min at reaction temperature (or as desired) to allow inhibitor/enzyme interaction.
12. Start assay by the addition of 50 μL H-Gly-Pro-pNA substrate (diluted in step 3) or 50 μL H-Gly-Pro-AMC substrate (diluted in step 4), which have been equilibrated to reaction temperature (e.g. 37°C). Enough of each substrate is provided to perform 96 assays at 100 μM for the chromogenic substrate or 5 μM for the fluorogenic substrate.
13. Read plate continuously, at $A_{405\text{nm}}$ for the pNA substrate, or Ex:380 nm/Em:460 nm for the AMC substrate, in a microplate reader. For example, record data at 1 min. intervals for a total of 10 to 60 min.
14. Perform data analysis (see below).

NOTE: Retain microplate for future use of unused wells.

TABLE 1. ASSAY MIXTURE EXAMPLES.

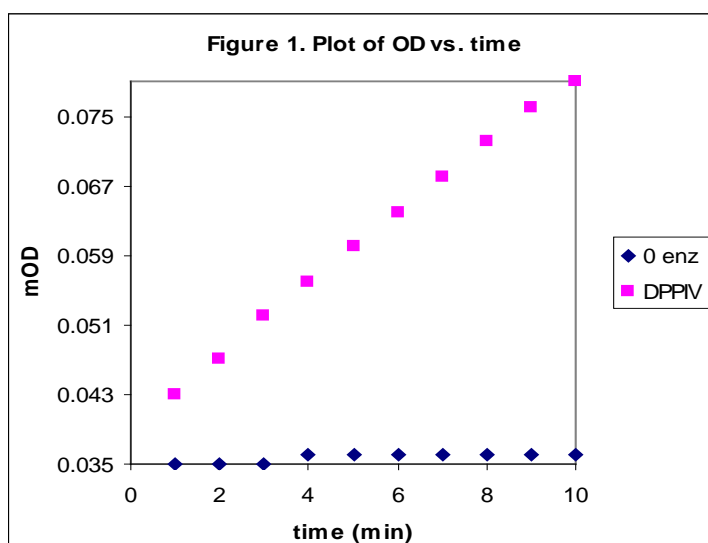
Sample	Assay Buffer	DPPIV (17.3 μU/μl)	Inhibitor	Substrate
Blank	50 μL	0	0	50 μL
Control	35 μL	15 μL	0	50 μL
Inhibitor [‡]	25 μL	15 μL	10 μL [‡]	50 μL
Test sample*	X μL	15 μL	Y μL*	50 μL

[‡]Refers to 10 μL of the diluted P32/98 prepared in step 2.

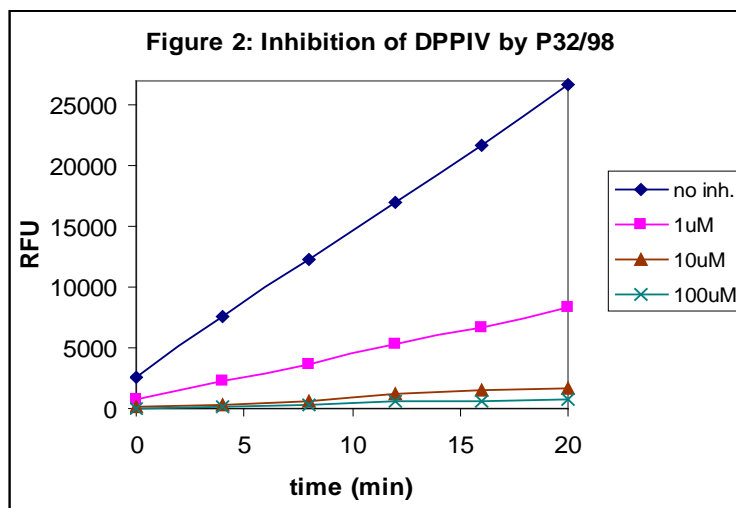
*Test sample is the experimental inhibitor. Dissolve/dilute experimental inhibitor into assay buffer and add to appropriate wells at desired volume “Y”. Adjust volume “X” to bring the total volume to 100 μL (X+Y=35 μL).

DATA ANALYSIS

- Plot data as A_{405nm} or Relative Fluorescence Units (RFU) versus time for each sample (see Figures 1 and 2).



Slope= $V=4.04E-03$ OD/min

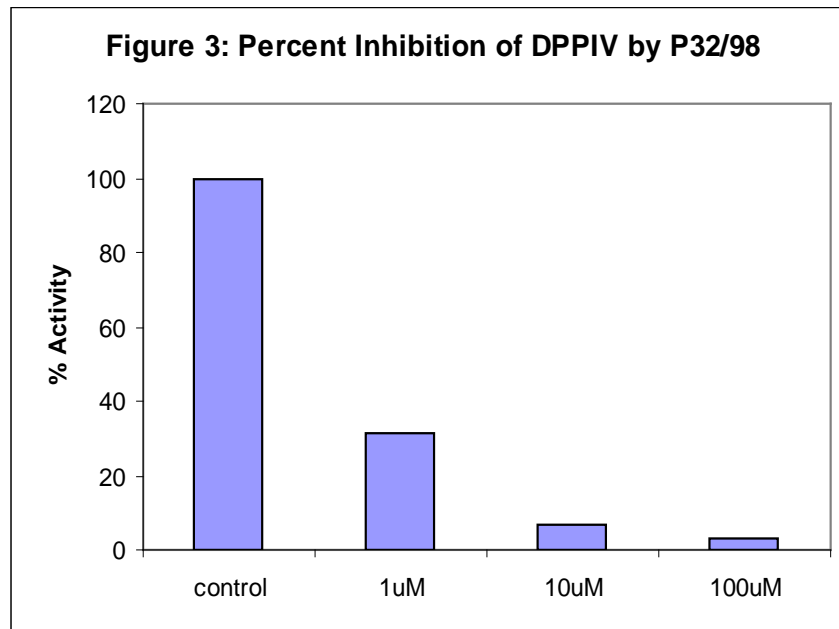


2. Determine the time range over which the reaction is linear. Typically, 0 to 15 min works well.
3. Obtain a “best fit” line for the data points and determine the slope.
4. Average the slopes of duplicate samples.
5. If the blank has a significant slope, subtract this number from the slopes for all samples.
6. To find % remaining activity in presence of inhibitor:

% activity remaining (with inhibitor)=

(slope of + inhibitor sample/control slope) x 100

See Figure 3.



7. To find the activity of the samples expressed as pmol substrate/min using chromogenic substrate:

Determine microplate reader conversion factor:

- a) Add 100 μL calibration standard (p-nitroaniline; 50 μM concentration) to 2 wells of the clear microplate. In Tris buffer the extinction coefficient for p-nitroaniline at 405 nm is $\sim 9700 \text{ M}^{-1}\text{cm}^{-1}$ (R. Lottenberg et al., 1983; *Biochim.Biophys.Acta* **742**:558). Typically, 100 μL of the 50 μM standard, in a $\frac{1}{2}$ volume well, produces an A_{405} of about 0.3. Alternatively, build a pNA standard curve and use the slope ($\mu\text{M}/\text{OD}$) as the conversion factor.
- b) Determine the average $A_{405\text{nm}}$ using 100 μL assay buffer as a blank.
- c) Calculate the conversion factor.

Conversion factor ($\mu\text{M}/\text{OD}$) = $50/\text{average } A_{405}$ from step b)

- d) Calculate the activity as pmol/min:

activity (pmol/min) = slope (OD/min) x conversion factor($\mu\text{M}/\text{OD}$) x assay vol (μL)

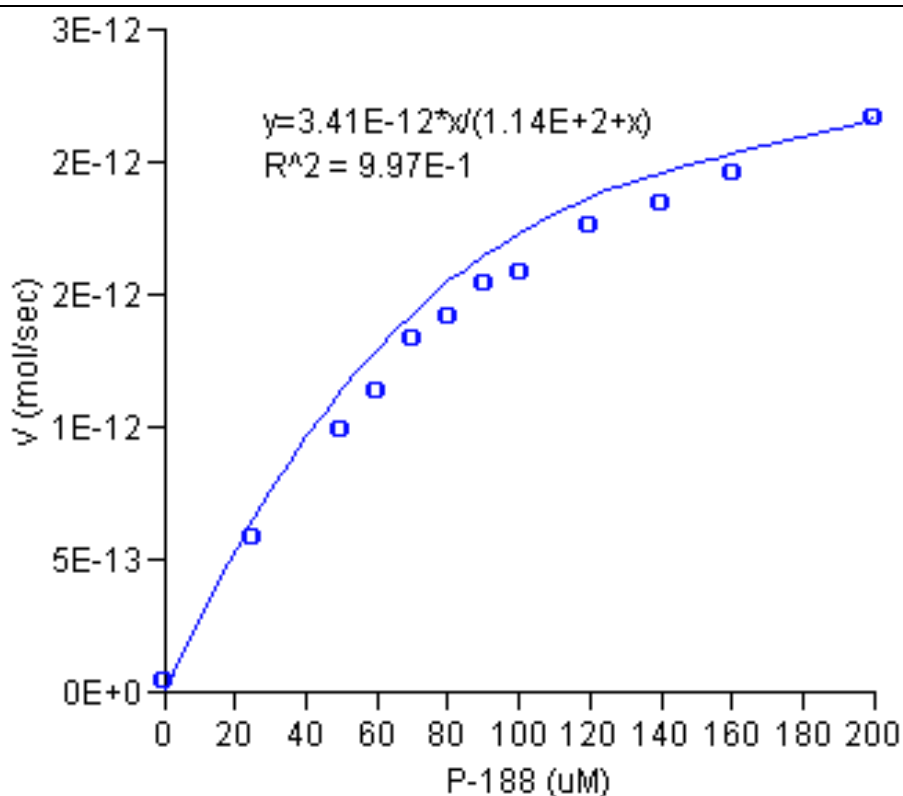
Example calculation for activity with colorimetric substrate:

conversion factor = $50 \mu\text{M}/0.294 \text{ OD} = 170 \mu\text{M}/\text{OD}$

activity of a control sample =

$4.04\text{E-}03 \text{ (OD/min)} \times 170(\mu\text{M}/\text{OD}) \times 100(\mu\text{l}) = 68 \text{ pmol/min}$

See Figure 4 for example of kinetic determination.



$$K_m=114 \mu\text{M}$$

$$V_{\text{max}}=3.41 \text{ pmol/sec}$$

$$k_{\text{cat}}/K_m=7.66 \times 10^5 \text{ M}^{-1}\text{s}^{-1}$$

Figure 4. Kinetics of H-Gly-Pro-pNA cleavage by DPPIV, 0.13 mU/well; 37°C. Rates were obtained from the slopes of the initial, linear portion of plots of A_{405} vs. time. Curve and kinetic parameters derived from a non-linear least squares fit to the Michaelis-Menten equation (Marquadt algorithm).

8. To find the activity of the samples expressed as pmol substrate/min using fluorogenic substrate:

a) Determine microplate reader conversion factor for AMC fluorophore. The exact AMC concentration range that will be useful for preparing a standard curve will vary depending on the fluorimeter model, the gain setting, and the exact excitation and emission wavelengths used. The AMC standard, as provided (30 μM), may yield off-scale readings in some cases. We recommend diluting some of the standard to a relatively low concentration with Assay Buffer (0.5 or 1.0 μM) and then measuring the fluorescence of 100 μl . The estimate of $\mu\text{M}/\text{RFU}$ obtained with this measurement, together with the observed range of values obtained in the enzyme assays, can then be used to plan an appropriate series of dilutions for a standard curve. The slope of the standard curve can then be used as the $\mu\text{M}/\text{RFU}$ conversion factor.

b) Calculate the activity as pmol/min:

$$\text{activity (pmol/min)} = \text{slope (RFU/sec)} \times 60\text{sec/min} \times \text{conversion factor}(\mu\text{M}/\text{RFU}) \times \text{assay vol } (\mu\text{L})$$

9. Inhibition kinetics:

For determining K_i , etc., the concentration of enzyme active sites is needed. Although the concentration (as measured by the Bradford method) of DPPIV in this kit can be obtained from Enzo Life Sciences Technical Services, it is more accurate to use active site titration with an inhibitor, such as P32/98 in this kit. Please see the following references:

a) *Active-site titration of peptidases*: C.G. Knight; *Methods Enzymol.* **248**, 85 (1995).

b) *L-trans-Epoxy succinyl-leucylamido(4-guanidino)butane (E-64) and its analogues as inhibitors of cysteine proteinases including cathepsins B, H and L*: A.J. Barrett et al.; *Biochem. J.* 201, 189 (1982).

NOTES

NOTES

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

www.enzolifesciences.com
Enabling Discovery in Life Science®

Global Headquarters

Enzo Life Sciences Inc.

10 Executive Blvd
Farmingdale, NY 11735
(p) 1-800-942-0430
(f) 1-631-694-7501
(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach
CH-4415 Lause / Switzerland
(p) +41/0 61 926 89 89
(f) +41/0 61 926 89 79
(e) info-ch@enzolifesciences.com

Please visit our website at www.enzolifesciences.com for additional contact information.