

**Product ID: K5100****Lot No.: 2110281**

Original Expiration Date: 19 July 2022

Extended Expiration Date*: 30 September 2024

Content	Concentration in Incubation
NADPH	0.675 mM

*: The expiration date was extended based on ongoing stability testing of the product.

NADPH is added to microsomal incubations to support reactions catalyzed by cytochrome P450, flavin-containing monooxygenases and certain other enzymes. RapidStart provides the required NADPH for these reactions by generating NADPH *in situ* through an enzymatic reaction catalyzed by glucose-6-phosphate.

Vial contents: 0.1 mL of 100 mM NADP
0.1 mL of 500 mM Glucose-6-phosphate
0.1 mL of 100 units/mL Glucose-6-phosphate dehydrogenase
In 20 mM potassium phosphate buffer, pH 7.4

Preparation procedures:

1. Determine the desired initial concentration using the dilution chart provided (page 2).
2. Add the appropriate amount of high purity water for desired concentration.
3. Vortex vial until all three pellets have fully dissolved.
4. NADPH has been generated and the working solution is ready for use.

Stability of activated NADPH regenerating system:

At 2-8°C: 48 hours

At room temperature: 12 hours

Do not refreeze.

The total number of incubations that can be performed is dependent on the assay design.

Suggested Use:**Store at -20°C or below***Intended for in vitro use only.*

These data were generated by and are the property of XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

DATASHEET PREPARED 26 MAY 2022

In a typical microsomal incubation, the recommended concentration of the NADPH-regenerating system consists of 1 mM NADP(H), 5 mM glucose-6-phosphate and 1 Unit/mL of glucose-6-phosphate dehydrogenase. These concentrations can be achieved by adding 0.70 mL of high purity water to the frozen components in the RapidStart vial, vortexing, and then adding the activated RapidStart solution to the incubation mixture in a 1:10 (v/v) dilution to initiate metabolic reactions.

RapidStart is designed as a flexible prepackaged NADPH regenerating system to accommodate various experimental designs. Depending on the amount of high purity water added to the system, RapidStart generates NADPH with a concentration of ~4-80 mM.

RapidStart Dilution Table

High Purity Water Added (mL)	NADP (mM)		G6P (mM)		G6PDH (units/mL)		Total Volume (mL)	Expected NADPH Concentration, As Prepared (mM)
	As Prepared	Final Incubation 1:10 (v/v)	As Prepared	Final Incubation 1:10 (v/v)	As Prepared	Final Incubation 1:10 (v/v)		
0.10	25.00	2.50	125.00	12.500	25.00	2.500	0.40	~17-22
0.20	20.00	2.00	100.00	10.000	20.00	2.000	0.50	~14-18
0.30	16.67	1.67	83.33	8.333	16.67	1.667	0.60	~10-13
0.40	14.29	1.43	71.43	7.143	14.29	1.429	0.70	~8.8-12.8
0.50	12.50	1.25	62.50	6.250	12.50	1.250	0.80	~10.0-11.2
0.60	11.11	1.11	55.56	5.556	11.11	1.111	0.90	~7.8-10
0.70*	10.00	1.00	50.00	5.000	10.00	1.000	1.00	~7.0-9.0
0.80	9.09	0.91	45.45	4.545	9.09	0.909	1.10	~6.4-8.2
0.90	8.33	0.83	41.67	4.167	8.33	0.833	1.20	~5.8-7.5
1.00	7.69	0.77	38.46	3.846	7.69	0.769	1.30	~5.4-6.9
1.10	7.14	0.71	35.71	3.571	7.14	0.714	1.40	~5.0-6.4
1.20	6.67	0.67	33.33	3.333	6.67	0.667	1.50	~4.7-6.0
1.30	6.25	0.63	31.25	3.125	6.25	0.625	1.60	~4.4-5.6
1.40	5.88	0.59	29.41	2.941	5.88	0.588	1.70	~4.1-5.3
1.50	5.56	0.56	27.78	2.778	5.56	0.556	1.80	~3.9-5.0
1.60	5.26	0.53	26.32	2.632	5.26	0.526	1.90	~3.7-4.7
1.70	5.00	0.50	25.00	2.500	5.00	0.500	2.00	~3.5-4.5

* Denotes XenoTech's recommended concentration.

NADPH Concentration Determination:

Although NADP and NADPH both absorb at 260 nm, only NADPH absorbs strongly at 340 nm. Consequently, the reduction of NADP to NADPH is associated with an increase in absorbance at ~340 nm. This increase in absorbance at ~340 nm can be used to determine the concentration of NADPH.

To verify the concentration of NADPH generated, a spectrophotometer is used to determine absorbance at ~340 nm (338-341nm is acceptable). Dispense 995 μ L of high purity water into a sample well and 1000 μ L into a reference well. Add 5 μ L of prepared RapidStart to the sample well and record the absorbance at 340 nm. The concentration of NADPH in an incubation can be determined using the following equation.

$$\frac{\text{Absorbance}}{\epsilon} \times \frac{\text{Volume of NADPH added to incubation mixture } (\mu\text{L})}{\text{Incubation volume } (\mu\text{L})} \times \frac{\text{Total volume in sample well } (\mu\text{L})}{\text{Volume of NADPH added to well } (\mu\text{L})} = \text{Concentration in incubation (mM)}$$

where ϵ = extinction coefficient of NADPH = 6.22 mM⁻¹cm⁻¹ (which indicates that a 1 mM solution of NADPH would have an absorbance value of 6.22 in a 1 cm pathlength).

These data were generated by and are the property of XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.