

# PRODUCT INFORMATION



## $\alpha$ -D-Galactose-1-phosphate (potassium salt hydrate)

Item No. 17863

**Formal Name:** 1-(dihydrogen phosphate)  $\alpha$ -D-galactopyranose, dipotassium salt, hydrate

**Synonym:** Gal-1-P

**MF:**  $C_6H_{11}O_9P \cdot 2K [XH_2O]$

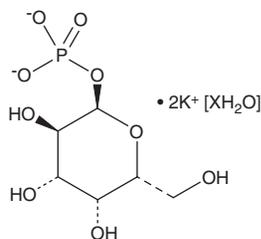
**FW:** 336.3

**Purity:**  $\geq 95\%$

**Supplied as:** A crystalline solid

**Storage:**  $-20^\circ\text{C}$

**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

$\alpha$ -D-Galactose-1-phosphate (potassium salt hydrate) is supplied as a crystalline solid. Aqueous solutions of  $\alpha$ -D-galactose-1-phosphate (potassium salt hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of  $\alpha$ -D-galactose-1-phosphate (potassium salt hydrate) in PBS (pH 7.2) is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

$\alpha$ -D-Galactose-1-phosphate is an intermediate in the interconversion of glucose and galactose. It is produced from galactose by galactokinase. Galactose-1-phosphate uridylyl transferase reversibly catalyzes the conversion of UDP-glucose (Item No. 15602) and galactose-1-phosphate to glucose-1-phosphate and UDP-galactose.

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/17/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM