

Xanthosine Kit

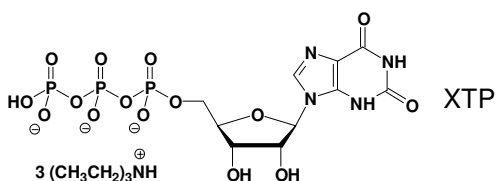
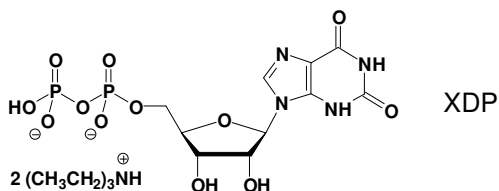
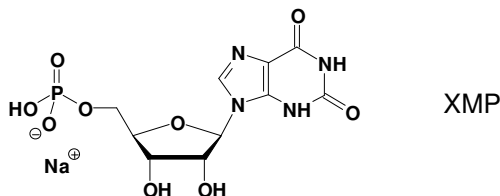
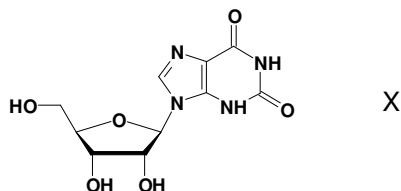
Cat.No.: NK-113

Kit Contents

Xanthosine Analogs	Cat. No.	Amount
X	N-1050	5 mg
XMP	NU-603	200 Units
XDP	NU-601	200 Units
XTP	NU-602	200 Units

1 Unit = 1 μ l of a 10 mM solution

Structures



Introduction

Xanthosines are used as G-nucleotide analogs in signal transduction research, molecular and cell biology.

Kit Description

The Xanthosine Kit contains a set of 4 typical Xanthosine analogs (Nucleoside, 5'-Mono, Di and Triphosphate).

Selected References:

Fulga *et al.* (2001) SR beta coordinates signal sequence release from SRP with ribosome binding to the translocon. *EMBO J.* **20** (9): 2338.

Legate *et al.* (2000) Nucleotide-dependent binding of the GTPase domain of the signal recognition particle receptor betasubunit to the alpha-subunit. *J. Biol. Chem.* **275** (35):27439.

Sakash *et al.* (2000) The use of nucleotide analogs to evaluate the mechanism of the heterotropic response of Escherichia coli aspartate transcarbamoylase. *Protein Sci.* **9** (1):53.

Muraoka *et al.* (1999) Effects of purinenucleotide analogues on microtubule assembly. *Cell Struct. Funct.* **24** (5):305.

Seifert *et al.* (1999) Effects of guanine, inosine, and xanthine nucleotides on beta(2)-adrenergic receptor/G(s) interactions: Evidence for multiple receptor conformations. *Mol. Pharmacol.* **56** (2):348.

Hwang *et al.* (1999) Structure-based identification of a novel NTPase from Methanococcus jannaschii. *Nat. Struct. Biol.* **6** (7):691.

Storage and Stability

Short term exposure (up to 1 week cumulative) to ambient temperature is possible. Long term storage is recommended at < -20°C. If properly stored, Jena Bioscience guarantees optimal performance of the compounds for 12 months after date of delivery.