

## UDP-Azide-Glucose

---

**Catalog No:** 55020

**Format:** 1 vial

**Quality Control:** 5-Hydroxymethylcytosine (5-hmC) is a DNA modification that results from the enzymatic conversion of 5-methylcytosine (5-mC) into 5-hydroxymethylcytosine by the TET family of cytosine oxygenases. One of the difficulties in studying 5-hmC is the fact that many of the traditional techniques employed to study DNA methylation do not distinguish between 5-mC and 5-hmC residues. In order to overcome some of the challenges with direct discrimination between 5-mC and 5-hmC, a T4 bacteriophage enzyme,  $\beta$ -glucosyltransferase, has been used to modify 5-hydroxymethylcytosine residues. The  $\beta$ -glucosyltransferase enzymes utilize a UDP-glucose (uridine diphosphoglucose) donor to attach a glucose moiety to 5-hmC which generates glucosyl-hydroxymethylcytosine. With the sugar attached to 5-hmC, the two types of DNA modifications can be discriminated using glucosyl-sensitive restriction enzymes or other techniques.

The UDP-Azide-Glucose contains a reactive group attached to the sugar. This enables the sugar to undergo further chemical modifications following treatment with the  $\beta$ -glucosyltransferase enzyme.

**Contents:** UDP-Azide-Glucose is provided at a concentration of 3 mM. Each vial contains 65  $\mu$ l.

**Storage and Guarantee:** Store at -20°C for up to 6 months. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.