

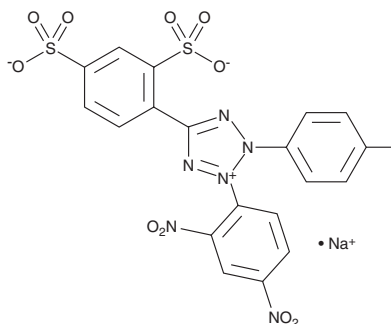
PRODUCT INFORMATION



WST-3

Item No. 30796

CAS Registry No.: 515111-36-1
Formal Name: 3-(2,4-dinitrophenyl)-5-(2,4-disulfophenyl)-2-(4-iodophenyl)-2H-tetrazolium, inner salt, monosodium salt
Synonym: Water-Soluble Tetrazolium 3
MF: $C_{19}H_{10}IN_6O_{10}S_2 \cdot Na$
FW: 696.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 233 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

WST-3 is supplied as a crystalline solid. A stock solution may be made by dissolving the WST-3 in water. The solubility of WST-3 in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

WST-3 is a water-soluble tetrazolium dye.¹ It is reduced by NADH in the presence of an electron mediator, such as 1-methoxy-5-methylphenazinium (1-methoxy PMS; Item No. 21258), to WST-3 formazan, which displays an absorption maxima of 433 nm.

Reference

1. Ishiyama, M., Sasamoto, K., Shiga, M., *et al.* Novel disulfonated tetrazolium salt that can be reduced to a water-soluble formazan and its application to the assay of lactate dehydrogenase. *Analyst* **120**, 113-116 (1995).

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, 11/09/2022

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