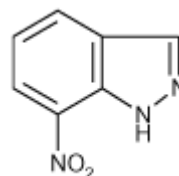


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 7-Nitroindazole (7-NI)**CATALOG #:** 00240**MOLECULAR INFORMATION:** C₇H₅N₃O₂
Mwt: 163.1
[2942-42-9]**PROPERTIES:****Color & Form** Yellow orange solid
Purity ≥ 98% by HPLC
Solubility Soluble in DMSO**STORAGE AND HANDLING:** Stored desiccated at 4°C upon receipt.**APPLICATION:** A selective inhibitor for the brain nitric oxide synthase (NOS).Ref: Moore, et al. *Methods in Enzymology* **268**, 393(1995); Faraci, et al. *Stroke* **26**(11), 2172(1995); Babbedge, et al. *Br. J. Pharmacol.* **110**(1), 225(1993).**TOXICITY:** Unknown**FIRST AID:** Potentially harmful; avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.