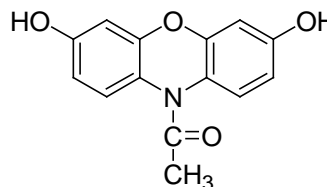


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 10-Acetyl-3,7-dihydroxyphenoxazine (*N*-Acetyl-3,7-dihydroxyphenoxazine)**CATALOG #:** 10061**MOLECULAR INFORMATION:** C₁₄H₁₁NO₄
Mwt: 257.24
[119171-73-2]**PROPERTIES:**

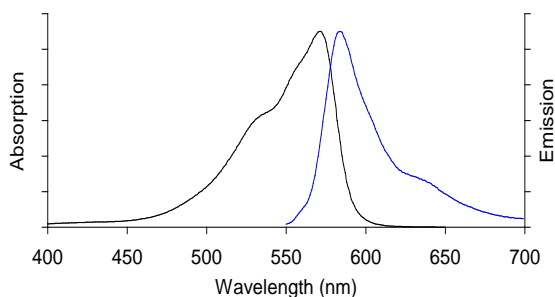
Color & Form	Off-white solid
Purity	≥ 98% by TLC
Solubility	Soluble in DMSO or DMF
Absorption/Emission	N/A
Extinction Coefficient	N/A

STORAGE AND HANDLING:

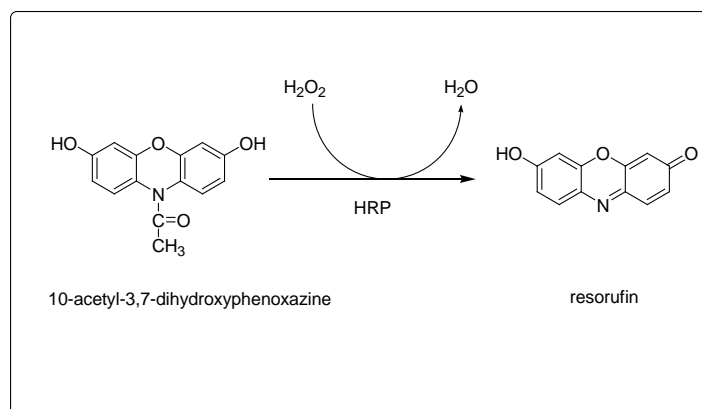
Store desiccated at -20 °C and protect from light.

APPLICATION:

10-Acetyl-3,7-dihydroxyphenoxazine, which has been marketed under the name Amplex Red™ reagent, is regarded as the best fluorogenic substrate for peroxidase because it is highly specific and stable. The substrate itself is nearly colorless and nonfluorescent until it is oxidized by H₂O₂ in the presence of horseradish peroxidase (HRP) to become the highly red fluorescent resorufin. High purity resorufin (#10062) is also available from Biotium to be used as a standard.

Ref.: 1) *J Immunol Methods* **202**, 133 (1997); 2) *Anal Biochem* **253**, 169(1997); 3) *Anal Biochem* **287**, 196 (2000)

Absorption and emission of resorufin in pH 8 buffer



Oxidation of 10-acetyl-3,7-dihydroxyphenoxazine forms highly fluorescent resorufin

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.*