

## DHE [Dihydroethidium]

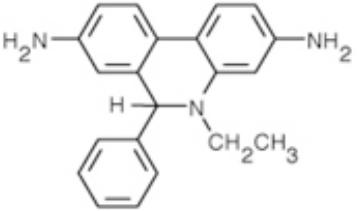
Catalog Number	Packaging Size
C260	25 mg

**Storage upon receipt:** -20°C, protected from light

### Introduction

Dihydroethidium (also called hydroethidium) is the chemically reduced form of the commonly used DNA dye ethidium bromide. The probe is useful to detect oxidative activities in viable cells, including respiratory burst in phagocytes. Dihydroethidium itself has blue fluorescence (Ex/Em: 355/420nm) in cells, while the oxidized form ethidium has red fluorescence (Ex/Em: 300/610 nm) upon DNA intercalation.

### Specifications

<b>Label:</b>	Ethidium	
<b>Ex/Em:</b>	300/610 nm	
<b>Detection Method:</b>	Fluorescent	
<b>Molecular Formula:</b>	C <sub>21</sub> H <sub>21</sub> N <sub>3</sub>	
<b>Molecular Weight:</b>	315.42	
<b>CAS Number:</b>	38483-26-0	
<b>Storage Conditions:</b>	-20°C, protected from light	
<b>Shipping Condition:</b>	Room Temperature	

### Applications

Probe for ROS

### References:

1. [G2019S leucine-rich repeat kinase 2 causes uncoupling protein-mediated mitochondrial depolarization.](#)  
 Papkovskaia TD, Chau KY, Inesta-Vaquera F, Papkovsky DB, Healy DG, Nishio K, Staddon J, Duchen MR, Hardy J, Schapira AH, Cooper JM,  
*Hum Mol Genet* (2012) 21:4201-4213
2. [Early safety assessment using cellular systems biology yields insights into mechanisms of action.](#)  
 Giuliano KA, Gough AH, Taylor DL, Vernetti LA, Johnston PA,  
*J Biomol Screen* (2010) 15:783-797
3. [Oxidant generation by single infected monocytes after short-term fluorescence labeling of a protozoan parasite.](#)  
 Chang HK, Thalhofer C, Duerkop BA, Mehling JS, Verma S, Gollob KJ, Almeida R, Wilson ME  
*Infect Immun* (2007) 75:1017-1024