

# PRODUCT DATA SHEET



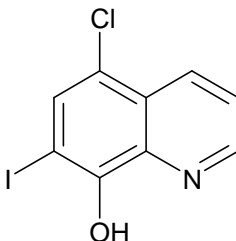
PRODUCT: Clioquinol  
5-Chloro-7-iodo-8-quinolinol

CAS NO.: 130-26-7

CATALOG NO.: N-158

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: Tan solid

MOLECULAR FORMULA:  $C_9H_5ClINO$

MOLECULAR WEIGHT: 305.5

PURITY: 98% (TLC: 5% Methanol/methylene chloride + 0.5%  $NH_4OH$ ;  $R_f = 0.25$ )

SOLUBILITY: Soluble in: DMSO (>25 mg/mL)

STORAGE: Store as supplied, at room temperature for up to 1 year. Store solutions at  $-20^\circ C$  for up to 3 months.

APPLICATION NOTES: A metal chelating ligand which displays protective effects in various neurodegenerative states. It has shown to modulate amyloid beta-peptide accumulation<sup>1</sup> via upregulation of metalloprotease activity<sup>2</sup>. When co-administered with vitamin B12 it totally rescues lead-induced impairments of synaptic plasticity in an animal model<sup>3</sup>. Clioquinol down-regulates mutant huntingtin expression *in vitro* and mitigates pathology in a Huntington's disease mouse model<sup>4</sup>.

- REFERENCES:
1. A. Caragounis *et al.* *Biochem. J.* 2007 **407** 435
  2. A.R. White *et al.* *J. Biol. Chem.* 2006 **281** 17670
  3. W.H. Chen *et al.* *Neuroscience* 2007 **147** 853
  4. T. Nguyen *et al.* *Proc. Natl. Acad. Sci. USA* 2005 **102** 11840

**The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.**