

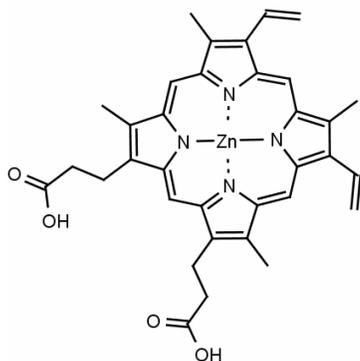


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Product Name: Zn (II) Protoporphyrin IX



Catalog Number	CAS Registry No.	MW (g/mol)	Chemical Formula	Solubility
F-H060	15442-64-5	626.03	C ₃₄ H ₃₂ N ₄ O ₄ Zn	Pyridine, DMF or DMSO

Storage: Zn (II) Protoporphyrin IX is stable for at least one year when stored as a solid, protected from moisture, at -20°C. Protect from light.

Field of Use: Zn (II) Protoporphyrin IX is an inhibitor of heme oxygenase¹ (the enzyme that catalyzes the conversion of heme to biliverdin in the heme degradation pathway) and guanylyl cyclase². Heme oxygenase has been implicated in tumor cell resistance to chemotherapy³, reduction of free radical formation⁴ and inflammation⁵ and associated with vascular repair⁵.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

References:

1. Maines, M. D. Zinc . protoporphyrin is a selective inhibitor of heme oxygenase activity in the neonatal rat. *Biochim Biophys Acta* **1981**, 673, 339-50.
2. Gupta, G.; Kim, J.; Yang, L.; Sturley, S. L.; Danziger, R. S. Expression and purification of soluble, active heterodimeric guanylyl cyclase from baculovirus. *Protein Expr Purif* **1997**, 10, 325-30.
3. Jozkowicz, A.; Was, H.; Dulak, J. Heme oxygenase-1 in tumors: is it a false friend? *Antioxid Redox Signal* **2007**, 9,2099-117.
4. Abraham, N. G.; Kappas, A. Heme oxygenase and the cardiovascular-renal system. *Free Radic Biol Med* **2005**, 39, 1-25.
5. Kim, D. H.; Burgess, A. P.; Li, M.; Tsenovoy, P. L.; Addabbo, F.; McClung, J. A.; Puri, N.; Abraham, N. G. Heme oxygenase-mediated increases in adiponectin decrease fat content and inflammatory cytokines, TNF and IL-6, in Zucker rats and reduce adipogenesis in human mesenchymal stem cells. *J Pharmacol Exp Ther* **2008**.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

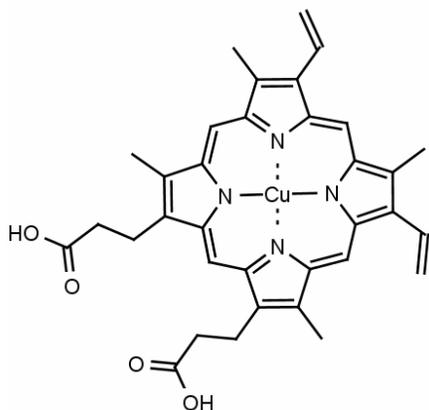


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Product Name: Cu (II) Protoporphyrin IX



Catalog Number	MW (g/mol)	Chemical Formula	Solubility
F-H070	624.19	C ₃₄ H ₃₂ CuN ₄ O ₄	Pyridine, DMF or DMSO

Storage: Cu (II) Protoporphyrin IX is stable for at least one year when stored as a solid, protected from moisture, at -20°C. Protect from light.

Field of Use: Cu (II) Protoporphyrin IX does not inhibit heme oxygenase (the enzyme that catalyzes the conversion of heme to biliverdin in the heme degradation pathway) and is used as a negative control for Zn (II) Protoporphyrin (an inhibitor of heme oxygenase). Heme oxygenase has been implicated in tumor cell resistance to chemotherapy¹, reduction of free radical formation² and inflammation³ and associated with vascular repair³.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

References:

1. Jozkowicz, A.; Was, H.; Dulak, J. Heme oxygenase-1 in tumors: is it a false friend? *Antioxid Redox Signal* **2007**, 9,2099-117.
2. Abraham, N. G.; Kappas, A. Heme oxygenase and the cardiovascular-renal system. *Free Radic Biol Med* **2005**, 39, 1-25.
3. Kim, D. H.; Burgess, A. P.; Li, M.; Tsenovoy, P. L.; Addabbo, F.; McClung, J. A.; Puri, N.; Abraham, N. G. Heme oxygenase-mediated increases in adiponectin decrease fat content and inflammatory cytokines, TNF and IL-6, in Zucker rats and reduce adipogenesis in human mesenchymal stem cells. *J Pharmacol Exp Ther* **2008**.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

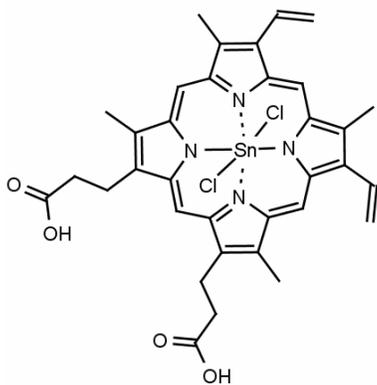


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Product Name: Sn (IV) Protoporphyrin IX Dichloride



Catalog Number	MW (g/mol)	Chemical Formula
F-H080	750.26	C ₃₄ H ₃₂ N ₄ O ₄ SnCl ₂

Storage: Sn (IV) Protoporphyrin IX is stable for at least one year when stored as a solid, protected from moisture, at -20°C. Protect from light.

Field of Use: Sn (IV) Protoporphyrin IX is an inhibitor heme oxygenase (the enzyme that catalyzes the conversion of heme to biliverdin in the heme degradation pathway) but has also been found to stimulate production of the heme oxygenase protein¹. Contrast the activity of Co (III) Protoporphyrin which has been found to have similar activities to Sn (IV) Protoporphyrin but with a greater enhancement of heme oxygenase synthesis activity such that heme oxygenase activity is actually increase when administered in vivo while in vitro administration inhibits heme oxygenase activity¹. Heme oxygenase has been implicated in tumor cell resistance to chemotherapy², reduction of free radical formation³ and inflammation⁴ and associated with vascular repair⁴.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

References:

1. Sardana, M. K.; Kappas, A. Dual control mechanism for heme oxygenase: tin(IV)-protoporphyrin potently inhibits enzyme activity while markedly increasing content of enzyme protein in liver. *Proc Natl Acad Sci U S A* **1987**, 84, 2464-8.
2. Jozkowicz, A.; Was, H.; Dulak, J. Heme oxygenase-1 in tumors: is it a false friend? *Antioxid Redox Signal* **2007**, 9,2099-117.
3. Abraham, N. G.; Kappas, A. Heme oxygenase and the cardiovascular-renal system. *Free Radic Biol Med* **2005**, 39, 1-25.
4. Kim, D. H.; Burgess, A. P.; Li, M.; Tsenovoy, P. L.; Addabbo, F.; McClung, J. A.; Puri, N.; Abraham, N. G. Heme oxygenase-mediated increases in adiponectin decrease fat content and inflammatory cytokines, TNF and IL-6, in Zucker rats and reduce adipogenesis in human mesenchymal stem cells. *J Pharmacol Exp Ther* **2008**.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

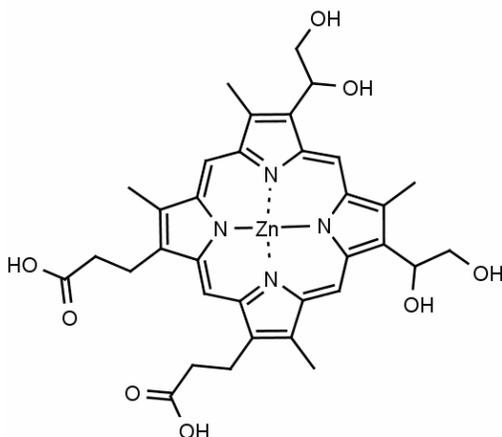


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Product Name: Zn (II) Deuteroporphyrin IX 2,4 bis ethyleneglycol



Catalog Number	MW (g/mol)	Chemical Formula	Solubility
F-H090	694.04	C ₃₄ H ₃₆ N ₄ O ₈ Zn	DMSO

Storage: Zn (II) Deuteroporphyrin IX 2,4 bis ethyleneglycol is stable for at least one year when stored as a solid, protected from moisture, at -20 °C. Protect from light.

Field of Use: Zn (II) Deuteroporphyrin IX 2,4 bis ethyleneglycol is an inhibitor of heme oxygenase (the enzyme that catalyzes the conversion of heme to biliverdin in the heme degradation pathway) at concentrations as low as 0.5 μM¹. Heme oxygenase has been implicated in tumor cell resistance to chemotherapy², reduction of free radical formation³ and inflammation⁴ and associated with vascular repair⁴.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

References:

1. Chernick, R. J.; Martasek, P.; Levere, R. D.; Margreiter, R.; Abraham, N. G. Sensitivity of human tissue heme oxygenase to a new synthetic metalloporphyrin. *Hepatology* **1989**, 10, 365-9.
2. Jozkowicz, A.; Was, H.; Dulak, J. Heme oxygenase-1 in tumors: is it a false friend? *Antioxid Redox Signal* **2007**, 9, 2099-117.
3. Abraham, N. G.; Kappas, A. Heme oxygenase and the cardiovascular-renal system. *Free Radic Biol Med* **2005**, 39, 1-25.
4. Kim, D. H.; Burgess, A. P.; Li, M.; Tsenovoy, P. L.; Addabbo, F.; McClung, J. A.; Puri, N.; Abraham, N. G. Heme oxygenase-mediated increases in adiponectin decrease fat content and inflammatory cytokines, TNF and IL-6, in Zucker rats and reduce adipogenesis in human mesenchymal stem cells. *J Pharmacol Exp Ther* **2008**.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.