



Dreiding stereomodels

Highly accurate, 3-dimensional, stainless steel molecular models for demonstrating stereochemical factors in conformational analysis and in the study of reaction mechanisms. Model units consist of push-together rods and tubes radiating from, and connecting to, a point that represents the atomic nucleus. Rods and tubes rotate to show conformational changes. Elements, angle strain, and bond types are identified by color-coded marks on the rods and tubes.

Set type	Total parts	Cat. No.
Simple	87	Z24,786-3
Normal	182	Z24,787-1
5-ring supplemental	51	Z24,789-8
3-ring supplemental	10	Z24,790-1
4-ring supplemental	30	Z24,792-8
7-ring supplemental	34	Z24,793-6
Peptide-A-Set	82	Z25,609-9
Porphyrin-Set	130	Z25,644-7
Nucleic-Acid-Set	393	Z25,645-5

Sets Contents

All models supplied in storage boxes.

Abbreviations:

Dobarom	Aromatic Double Bond	plan	Planar
Dobcon	Conjugated Double Bond	frot	Free Rotating
Dob (Dobisol)	Isolated Double Bond	lin	Linear
endo	Endocyclic	*	unit in which one or more rods are substituted by tubes
exo	Exocyclic		

Simple-Set (15 58 000-3) Z24,786-3

- 2 ACETYLENE
- 5 BIOXYGEN
- 5 CARBONYL
- 9 DOBAROM
- 5 DOBCON
- 5 DOBISOL
- 50 TETRACARBON
- 5 TRINITROGEN
- 1 ANGSTROMSCALE

Normal-Set (15 58 010-1) Z24,787-1

- 2 ACETYLENE
- 2 AMIDE
- 2 BENZENE
- 5 BIOXYGEN
- 1 BISULFUR
- 5 CARBONYL
- 9 DOBAROM
- 5 DOBCON
- 5 DOBISOL
- 2 IMINE
- 1 *IMINE
- 1 IMONIUM

- 1 MONOBROMO
- 1 *MONOBROMO
- 1 MONOCHLORO
- 1 *MONOCHLORO
- 1 MONOFLUORO
- 1 *MONOFLUORO
- 1 MONOIODO
- 1 *MONOIODO
- 1 MONOXYGEN
- 1 *MONOXYGEN
- 1 NITRILE
- 1 *NITRILE
- 50 TETRACARBON
- 2 TETRANITROGEN

- 1 TETRAPHOSPHOROUS
- 1 TRICARBONPLAN
- 1 *TRICARBONPLAN
- 5 TRINITROGEN
- 2 TRINITROGENPLAN
- 1 TRIPHOSPHOROUS
- 1 ANGSTROMSCALE
- 12 C-EXTENSION
- 12 H-EXTENSION
- 12 *H-EXTENSION
- 20 ROTORESISTOR
- 5 VAN DER WAALS H-RADIUM
- 5 *VAN DER WAALS H-RADIUS

5-Ring-Supplement-Set (15 58 020-8) Z24,789-8

- 1 AMIDE5ENDO
- 1 *AMIDE5ENDO
- 1 BISULFUR5
- 3 CARBONYL5
- 4 DOBAROM5ENDO
- 4 DOBCON5ENDO
- 1 DOBCON5ENDO5EXO
- 1 DOBCON5EXO
- 1 DOBCON5EXO6ENDO
- 5 DOB5ENDO

- 3 DOB5EXO
- 1 IMINE5CEXO
- 1 IMINE5ENDO
- 1 *IMINE5ENDO
- 1 IMONIUM5ENDO
- 1 *IMONIUM5ENDO
- 1 IMONIUM5NEXO
- 15 TETRACARBON5
- 3 *TETRACARBON5
- 1 TRINITROGEN5PLAN
- 1 *TRINITROGEN5PLAN

3-Ring-Supplement-Set (15 58 030-5) Z24,790-1

- 4 CYCLOPROPANE
- 3 ETHYLENE OXIDE
- 3 ETHYLENEIMINE

4-Ring-Supplement-Set (15 58-040-2) Z24,792-8

- 1 AMIDE4ENDO
- 1 BIOXYGEN4
- 2 CARBONYL4
- 3 CARBONYL4
- 3 DOBAROM4ENDO
- 3 DOB4ENDO
- 2 DOB4EXO
- 12 TETRACARBON4
- 5 *TETRACARBON4
- 1 TRINITROGEN4

**7-Ring-Supplement-Set
(15 58 050-0)**

Z24,793-6

- 1 BIOXYGEN7
- 1 CARBONYL7
- 20 TETRACARBON7
- 5 *TETRACARBON7
- 5 TRICARBONCON7PLAN
- 2 *TRICARBONCON7PLAN

**Peptide-A-Set (15 58 060-7)
Z25,609-9**

- 60 AMIDE
- 6 BENZENE
- 10 BIOXYGEN
- 6 BISULFUR
- 5 CARBONYL
- 6 DOBAROM
- 3 DOBAROM5ENDO
- 3 DOBCON5ENDO
- 3 DOB5ENDO
- 1 GUANIDINIUM
- 3 IMINE5ENDO
- 60 INTERHYDROGEN
(Hydrogen Bond)
- 5 MONOXYGEN
- 1 PROLINE (TRANS)
- 150 TETRACARBON
- 5 TETRANITROGEN
- 3 TRINITROGEN5PLAN
- 1 ANGSTROMSCALE

**Porphyrin-Set (15 58 070-4)
Z25,644-7**

- 2 AMIDE
- 2 *AMIDE
- 6 BINITROGEN5
- 5 BIOXYGEN
- 5 CARBONYL
- 12 DOBCON5ENDO
- 3 DOBCON5ENDO5EXO
- 6 DOBCON5EXO
- 5 DOBISOL
- 2 HEXACOBALT
- 2 HEXAIRON
- 3 IMONIUM5ENDO
- 3 *IMONIUM5ENDO
- 20 TETRACARBON
- 2 TETRANICKELPLAN
- 12 TRICARBONCON5PLAN
- 12 *TRICARBONCON5PLAN
- 12 TRICARBONCON7PLAN
- 3 *TRICARBONCON7PLAN
- 6 TRINITROGEN5PLAN
- 6 *TRINITROGEN5PLAN
- 1 ANGSTROMSCALE

**Nucleic-Acid-Set (15 58 080-1)
Z25,645-5**

(example: DNA with 5 Thymine, 5 Adenine, 6 Cytosine, 6 Guanine)

- 5 AMIDE
- 44 BIOXYGEN
- 44 BIOXYGEN7
- 17 CARBONYL
- 11 DOBAROM
- 11 DOBAROM5ENDO
- 6 IMINE
- 5 *IMINE
- 6 IMINE5ENDO
- 5 *IMINE5ENDO
- 11 IMONIUM
- 20 MONOXYGEN
- 113 TETRACARBON
- 20 TETRAPHOSPHOROUS
- 17 TRINITROGENPLAN
- 17 *TRINITROGENPLAN
- 11 *TRINITROGEN5PLAN
- 11 BASE-PAIR-SUPPORT
- 1 DNA-BASE WITH ROD
- 17 DNA-H-BOND
- 1 ANGSTROMSCALE



ALDRICH®

*Organics and Inorganics for
Chemical Synthesis*

1001 West Saint Paul Ave., Milwaukee, WI 53233

Telephone 414-273-3850 • 800-231-8327

Fax 414-273-4979 • 800-962-9591

Internet <http://www.sial.com/aldrich>

E-mail aldrich@sial.com

© 1997 Sigma-Aldrich Co. Printed in the USA. Aldrich is a member of the Sigma-Aldrich Family. Aldrich warrants that its products conform to the information contained in this and other Aldrich publications. Purchaser must determine the suitability of the product for its particular use. See reverse side of invoice or packing slip for additional terms and conditions of sale.