

AL-231

Technical Bulletin

Properties of Elastomers Chart

PROPERTIES:	Ethylene Propylene	Fluorocarbon	Natural Rubber	Neoprene	Nitrile	Silicone	Styrene Butadiene	Urethane Rubber
Tear Resistance	Good	Fair	Good	Good	Fair	Poor	Fair	Excellent
Abrasive Resistance	Good	Good	Excellent	Excellent	Good	Poor	Good	Outstanding
Compression Set	Good	Very Good	Good	Fair to Good	Good	Fair	Excellent	Good
Permability to Gases	Fair to Poor	Very Low	Fair	Low	Fair	Fair	Fair	Fair
ACID RESISTANCE:								
Dilute	Excellent	Excellent	Fair to Good	Very Good	Good	Fair	Fair to Good	Poor
Concentrated	Good	Excellent	Fair to Good	Good	Good	Fair	Fair to Good	Poor
SOLVENT RESISTANCE:								
Aliphatic Hydrocarbons	Poor	Excellent	Poor	Good	Excellent	Poor	Poor	Excellent
Aromatic Hydrocarbons	Poor	Excellent	Poor	Fair	Good	Poor	Poor	Fair to Good
Oxygenated (ketones, etc.)	Good	Poor	Good	Poor	Poor	Fair	Poor	Poor
RESISTANCE:								
Oil and Gasoline	Poor	Excellent	Poor	Good	Excellent	Fair	Poor	Excellent
Water Absorption	Excellent	Very Good	Very Good	Good	Fair to Good	Good	Very Good	Good @ 70F Poor @ 212 F
Ozone	Excellent	Excellent	Fair	Excellent	Fair	Excellent	Fair	Outstanding
Sunlight	Very Good	Very Good	Poor	Very Good	Poor	Excellent	Fair	Good
Temp. Range °C	-50 to 150	-20 to 200	-55 to 90	-40 to 100	-40 to 100	-60 to 200	-50 to 100	-25 to 100
Major Attributes	General Purpose	Heat Resistant, Excellent Solvent & Chemical Resistance	Resilience	Oil & Gas Resistance, Weatherability	Oil Resistance	Heat Resistance	General Purpose	Abrasion Resistance, Load Bearing Characteristics