

AL-224

Technical Bulletin

Description of Soda Lime Glass Used in Corning Labware

Soda Lime Glass
Code No. 0215
Used for Flint Glass

Standards

Glass commonly referred to as
Type II Soda Lime Glass per
A.S.T.M.-438 federal spec DD-G-54Ib

Chemical Composition

Composition	(percent approx.)
SiO ₂	73%
Na ₂ O	14%
CaO	7%
MgO	4%
Al ₂ O ₃	2%

Physical Properties

Coeff. of Exp.	89 x 10 ⁻⁷ cm/cm/°C
Strain Point	511°C
Anneal Point	545°C
Softening Point	724°C
Density	2.40 g/cm ³
Young's Mod.	Not available
Refract. Index	1.515 @ Sodium D Line

Transmittance/Wavelength For 0215 Soda Lime Glass - Approximately 1mm Thick

Applications

General purpose glass used in the container and plate glass industry.

Products

Microscope Slides and Pasteur Pipets.

Warnings

1. Do not use hydrofluoric or hot phosphoric acid on glass.
-