

AL-226

Technical Bulletin

Description of Zinc Titania Glass Used in Corning Labware

Zinc Titania Glass
Code No. 0211
Used for Cover Glass

Standards
Proprietary Corning composition.
No current federal or ASTM standards on this glass.

Chemical Composition

Composition	(percent approx.)
SiO ₂	64%
B ₂ O ₃	9%
ZnO	7%
K ₂ O	7%
Na ₂ O	7%
TiO ₂	3%
A ₁₂ O ₃	3%

Physical Properties

Coeff. of Exp.	74 x 10 ⁻⁷ cm/cm/°C
Strain Point	508°C
Anneal Point	550°C
Softening Point	720°C
Density	2.53 g/cm ³
Young's Mod.	7.6 x 10 ³ Kg/mm ²
Refract. Index	1.523 @ Sodium D Line
Surface Coating	Proprietary

Transmittance/Wavelength For 0211 Zinc Titania Glass - Approximately 1mm Thick

Applications

Designed specifically as a cover glass for attachment to microscope slides to preserve and protect samples on the slides. Also ideal for making blood smears suitable for microscopic examination.

Products

Cover glass and Micro-Sheet.

Warnings

1. Do not use in heat applications.