

# Calculating Nucleic Acid or Protein Concentration Using the GloMax<sup>®</sup>-Multi+ Microplate...

Calculating Nucleic Acid or Protein Concentration Using the GloMax<sup>®</sup>-Multi+ Microplate Instrument

---

## ABSTRACT

*Direct measurements of nucleic acid samples at OD<sub>260</sub> or protein samples at OD<sub>280</sub> can be converted to concentration using the Beer-Lambert law which relates absorbance to concentration using the pathlength of the measurement and an extinction coefficient. This application note discusses the relationship between these two variables and provides protocols quantifying proteins and nucleic acids.*



## Download Article

(229 KB)

---