

Gender Differences in Protein Kinase Expression

Mouse
Submandibular
Gland

