

Gender Differences in Protein Kinase Expression

Mouse Brain

