

Quanterix Board Diversity Matrix as of October 2, 2024

Quanterix Board Diversity Matrix as of October 2, 2024				
Total Number of Directors	9			
	Female	Male	Non-Binary	Did Not Disclose Gender
Part I: Gender Identity				
Directors	3	5	-	1
Part II: Demographic Background				
African American or Black	-	-	-	-
Alaskan Native or Native American	-	-	-	-
Asian	-	-	-	-
Hispanic or Latinx	-	-	-	-
Native Hawaiian or Pacific Islander	-	-	-	-
White	3	5	-	-
Two or More Races or Ethnicities	-	-	-	-
LGBTQ+				
Did Not Disclose Demographic Background	1			