

| Plate    | Position | Catalog No. | Mature miRNA Accession No. | Mature miRNA ID | miRNAs also Amplified |
|----------|----------|-------------|----------------------------|-----------------|-----------------------|
| QM063_01 | A01      | HmiRQP0098  | MIMAT0000444               | hsa-mir-126     |                       |
| QM063_01 | A02      | HmiRQP0099  | MIMAT0000445               | hsa-mir-126-3p  |                       |
| QM063_01 | A03      | HmiRQP0284  | MIMAT0000226               | hsa-miR-196a-5p |                       |
| QM063_01 | A04      | HmiRQP0285  | MIMAT0004562               | hsa-miR-196a-3p |                       |
| QM063_01 | A05      | HmiRQP0078  | MIMAT0005898               | hsa-miR-1246    |                       |
| QM063_01 | A06      | HmiRQP0378  | MIMAT0000688               | hsa-mir-301a    |                       |
| QM063_01 | A07      | HmiRQP0378  | MIMAT0000688               | hsa-mir-301a-3p |                       |
| QM063_01 | A08      | HmiRQP2236  | MIMAT0019704               | hsa-miR-4644    |                       |
| QM063_01 | A09      | HmiRQP1554  | MIMAT0016858               | hsa-miR-4306    |                       |
| QM063_01 | A10      | HmiRQP0195  | MIMAT0004608               | hsa-mir-146a    |                       |
| QM063_01 | A11      | HmiRQP0220  | MIMAT0004658               | hsa-mir-155     |                       |
| QM063_01 | A12      | HmiRQP0033  | MIMAT0004556               | hsa-mir-10b     |                       |
| QM063_01 | B01      | HmiRQP0315  | MIMAT0004494               | hsa-mir-21      |                       |
| QM063_01 | B02      | HmiRQP0509  | MIMAT0001631               | hsa-miR-451a    |                       |
| QM063_01 | B03      | HmiRQP2171  | MIMAT0019064               | hsa-miR-4525    |                       |
| QM063_01 | B04      | HmiRQP0316  | MIMAT0000076               | hsa-miR-21-5p   |                       |
| QM063_01 | B05      | HmiRQP0315  | MIMAT0004494               | hsa-miR-21-3p   |                       |
| QM063_01 | C01      | NC          | NC                         | NC              |                       |
| QM063_01 | C02      | NC          | NC                         | NC              |                       |
| QM063_01 | C03      | HKG1        | HKG1                       | HKG1            |                       |
| QM063_01 | C04      | HKG2        | HKG2                       | HKG2            |                       |
| QM063_01 | C05      | HKG3        | HKG3                       | HKG3            |                       |
| QM063_01 | C06      | HKG4        | HKG4                       | HKG4            |                       |
| QM063_01 | C07      | HKG5        | HKG5                       | HKG5            |                       |
| QM063_01 | C08      | HKG6        | HKG6                       | HKG6            |                       |
| QM063_01 | C09      | RT          | RT                         | RT              |                       |
| QM063_01 | C10      | RT          | RT                         | RT              |                       |
| QM063_01 | C11      | PCR         | PCR                        | PCR             |                       |
| QM063_01 | C12      | PCR         | PCR                        | PCR             |                       |