#### FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

	9	,		

l	OMB APPROVAL					
	OMB Number:	3235-0287				
l	Estimated average bu	rden				

0.5

hours per response:

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).

### STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(b) of the Investment Company Act of 1940

- D(-) t l								
g Person(s) to Issuer								
10% Owner								
Other (specify below)								
SVP, Commercial & Accelerator								
Filing (Check Applicable								
e Reporting Person								
re than One Reporting								
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned								
6. Ownership 7. Nature								
Form: Direct of Indirect (D) or Indirect Beneficial								
(I) (Instr. 4) Ownership (Instr. 4)								
D								
of 10. 11. Nature of Indirect								
Form: Beneficial								
Direct (D) Ownership or Indirect (Instr. 4)								
(I) (Instr. 4)								
(s)								
1 1								
1 1								

#### **Explanation of Responses:**

- 1. The sale of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 148 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$34.50 to \$34.8384. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.
- 3. Includes 12,281 restricted stock units.

# Remarks:

/s/ Brian Keane, Attorney-infact 07/03/2019

\*\* Signature of Reporting Person Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

- \* If the form is filed by more than one reporting person, see Instruction 4 (b)(v).
- \*\* Intentional misstatements or omissions of facts constitute Federal Criminal Violations See 18 U.S.C. 1001 and 15 U.S.C. 78ff(a).

Note: File three copies of this Form, one of which must be manually signed. If space is insufficient, see Instruction 6 for procedure.

Persons who respond to the collection of information contained in this form are not required to respond unless the form displays a currently valid OMB Number.