Instruction 1(b).

FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
-------------	------	-------

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL						
OMB Number: 3235-0287						
Estimated average burden						
hours per response:	0.5					

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

(Check all applicable) Director 10% Owner		Symbol		ker or Tr							f Reporting Person*	nd Address of David C.																			
A below) below)	3. Date of Earliest Transaction (Month/Day/Year)									Middle) N,	rst) (I CORPORATIOI TURNPIKE																				
onth/Day/Year) 6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person	Line) X Form filed by One Reporting Porm filed by More than One R					4. If A		1821 Zip)	A 0	ICA MA	(Street) BILLERI (City)																				
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																															
Securities Acquired (A) or posed Of (D) (Instr. 3, 4 and Posed Of (D) or Indirect (D) or Indirect (D) (Instr. 4) (I) (Instr. 4)				3. Transa Code (8)	Execution Date,			2. Transaction Date (Month/Day/Year		tr. 3)	Security (Inst	1. Title of S																			
(A) or Transaction(s)	(A) or PI	Amount	v	Code																											
,667 ⁽¹⁾ D \$43.04 42,844 ⁽²⁾ D	1,667 ⁽¹⁾ D \$		1,667 ⁽¹⁾ D \$		1,667 ⁽¹⁾ D					2020	10/16/	Common Stock																			
d of, or Beneficially Owned vertible securities)											Tal																				
Amount of Securities Underlying Derivative Security (Instr. 5) Derivative Security (Instr. 3 and 4) Derivative Security (Instr. 5) Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)	te Amount of Securities Underlying Derivative Security (Ins 3 and 4) Expiration Amount of Securities Security (Ins 3 and 4)		Expiration Dat (Month/Day/Ya		Expiration Date (Month/Day/Year		Expiration Date (Month/Day/Year)		Expiration Date (Month/Day/Year)		(Month/Day/Year)		Expiration Date (Month/Day/Year)		Expiration Date (Month/Day/Year)		Expiration Date (Month/Day/Year)		Expiration Date (Month/Day/Year)		Expiration Date Month/Day/Year) Solution Do Se 3 3		rities ired r osed)	5. Nu of Deriv Secu Acqu (A) or Dispo of (D) (Instr and 5		4. Transa Code (8)	emed ion Date, //Day/Year)	if any	3. Transaction Date (Month/Day/Year)	2. Conversion or Exercise Price of Derivative Security	1. Title of Derivative Security (Instr. 3)
Count (A) or (D) Price Transaction(s) (Instr. 3 and 4)	D \$ or Benefic le securiti 7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4) Amou or Numb	1,667 ⁽¹⁾ Dsed of, oonvertible isable and te ear)	Dispo ons, o e Exerc tion Da n/Day/Y	S sired, I optio 6. Date Expirat (Month	mber rative rities ired r osed	5. Nu of Deriv Secu Acqu (A) or Dispo of (D)	alls, v	ive Se Its, Ca 4. Transa Code (Derivat (e.g., pu	3A. Dee Execut if any	3. Transaction Date	2. Conversion or Exercise Price of Derivative	1. Title of Derivative Security																		

Explanation of Responses:

- 1. The sale of shares reported on this Form 4 was made pursuant to a Rule 10b5-1 trading plan adopted on May 15, 2020.
- 2. Includes 12,082 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

** Signature of Reporting Person

Date

10/20/2020

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.